

PDS-1000/He Biolistic Particle Delivery System

Instruction Manual

Catalog Number #1652257
#1652258

For technical service call your local Bio-Rad Office.
In the U.S. call 1-800-4BIORAD (1-800-424-6723).



BIO-RAD

Warranty and Regulatory Notices

Warranty Statement

The PDS-1000/He System and associated accessories are covered by a standard Bio-Rad warranty. Contact your local Bio-Rad Laboratories office for details of the warranty.

Regulatory Notices

Important: This Bio-Rad instrument is designed and certified to meet EN/IEC 61010-1:2010 (3rd Edition), EN61326-1:2013 Class A, EN50581:2012 requirements, which are internationally accepted electrical safety standards. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will void the manufacturer's warranty, void the regulatory certifications, and create a potential safety hazard.

Note: This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radiofrequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference, in which case the user will be required to correct the interference at his own expense.

Table of Contents

Section 1	Introduction to Particle Delivery	1
	Particle Delivery Technology	1
	Overview of PDS-1000/He Particle Delivery System	1
	Important Safety Information	2
	Requirements for System Operation	3
Section 2	Product Description.	5
	Packing List	5
	Identification of Unit Controls and Components	9
Section 3	Installation	11
	Connecting the PDS-1000/He System to a Helium Source	11
	Connecting the PDS-1000/He System to a Vacuum Source	13
	Power Cord/Voltage Regulator	14
Section 4	Operation of the PDS-1000/He Instrument	15
	Preparation of System Components Prior to Bombardment	15
	Performing a Bombardment	20
	Removal of Residual Helium Pressure – Shutdown	28
Section 5	Selection and Adjustment of System Bombardment Parameters	28
	Overview — Matrix of Variables	28
	Vacuum Level in Bombardment Chamber	29
	Helium Pressure/Rupture Disk Selection	29
	Solenoid Valve Adjustment	29
	Vacuum Flow Rate Control Valves	29
	Distance Between Rupture Disk and Macrocarrier	29
	Distance Between Macrocarrier and Stopping Screen	30
	Distance Between Stopping Screen and Target Shelf (Microcarrier Flight Distance)	30
	Microcarrier Selection	31
	Preparation of Biological Material for Bombardment	31
Section 6	Troubleshooting.	34
	Rupture Disk Bursts at Incorrect Pressure.	34
	Stopping Screen Forced through Screen Support Ring	34
	Excessive Gas Usage	35
	Chamber Will Not Hold Vacuum	35
	Sample Damage from Gas Pressure Wave	35
	Unit Will Not Pressurize Gas Acceleration Tube.	35
Section 7	Product Information.	36
	Biolistic Particle Delivery System.	36
	Spare Parts for Helium Biolistic System Assembly (PDS-1000/He)	36
Section 8	Appendices	37
	Cleaning the PDS-1000/He Device	37
	Metal Case Version	38
	Specifications	41
	Performing a Bombardment — Quick Guide Tear-Out	42
Section 9	References	43

Section 1

Introduction to Particle Delivery

For additional PDS-1000/He training and overview information, including instrument setup, consumable preparation, and instrument operation and shutdown, go to bio-rad.com/PDS1000/training.

Particle Delivery Technology

Biolistic particle delivery is a method of transformation that uses helium pressure to introduce DNA-coated microcarriers into cells. Microprojectile bombardment can transform such diverse targets as bacterial, fungal, insect, plant, and animal cells and intracellular organelles. Particle delivery is a convenient method for transforming intact cells in culture since minimal pre- or postbombardment manipulation is necessary. In addition, this technique is much easier and faster to perform than the tedious task of micro-injection. Both stable and transient transformations are possible with the biolistic particle delivery system.

Overview of PDS-1000/He Particle Delivery System

The Biolistic System

The biolistic PDS-1000/He Instrument uses pressurized helium to accelerate subcellular sized microprojectiles coated with DNA (or other biological material) over a range of velocities necessary to optimally transform many different cell types. The system consists of the bombardment chamber (main unit), connective tubing for attachment to a vacuum source, and all components necessary for attachment and delivery of high-pressure helium to the main unit (helium regulator, solenoid valve, and connective tubing).

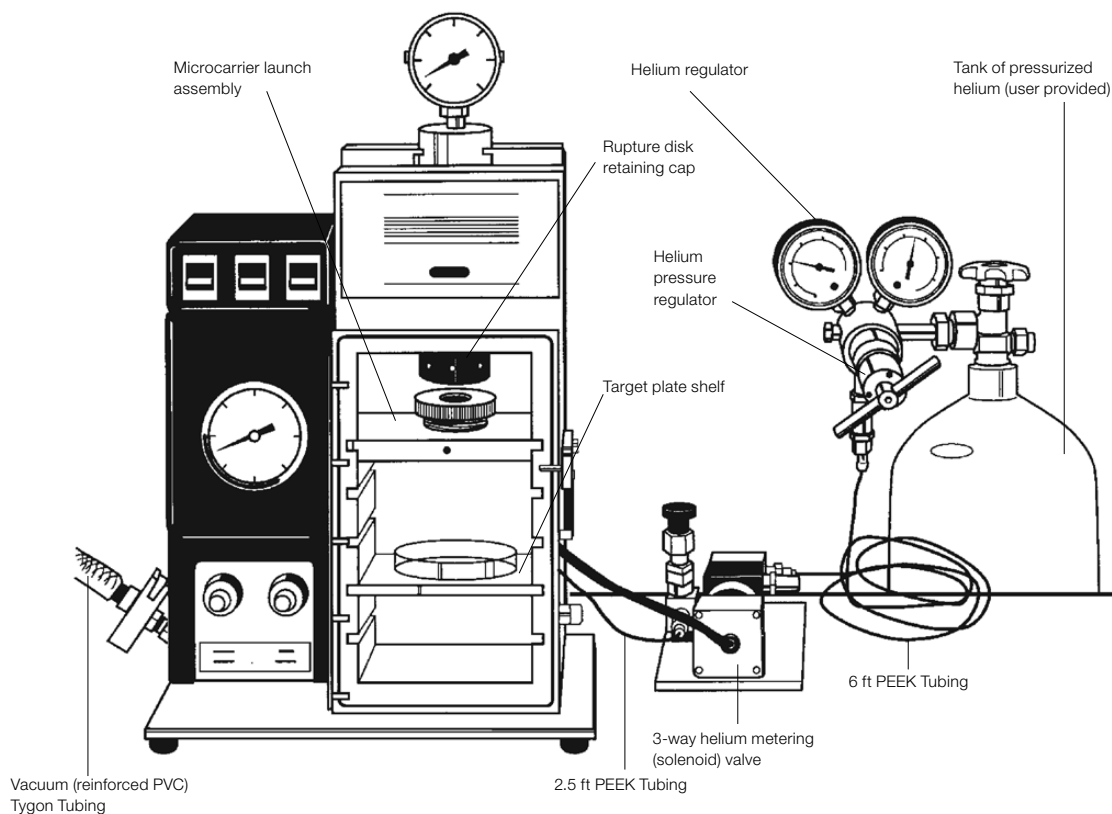


Fig. 1.1. Unit components, front view.

The Biolistic Process

The biolistic PDS-1000/He System uses high-pressure helium, released by a rupture disk, and partial vacuum to propel a macrocarrier sheet loaded with millions of microscopic gold microcarriers toward target cells at high velocity. The microcarriers are coated with DNA or other biological material for transformation. The macrocarrier is halted after a short distance by a stopping screen. The DNA-coated microcarriers continue traveling toward the target to penetrate and transform the cells.

The launch velocity of microcarriers for each bombardment depends on the helium pressure (rupture disk selection), the amount of vacuum in the bombardment chamber, the distance from the rupture disk to the macrocarrier (A), the macrocarrier travel distance to the stopping screen (B), and the distance between the stopping screen and target cells (C).

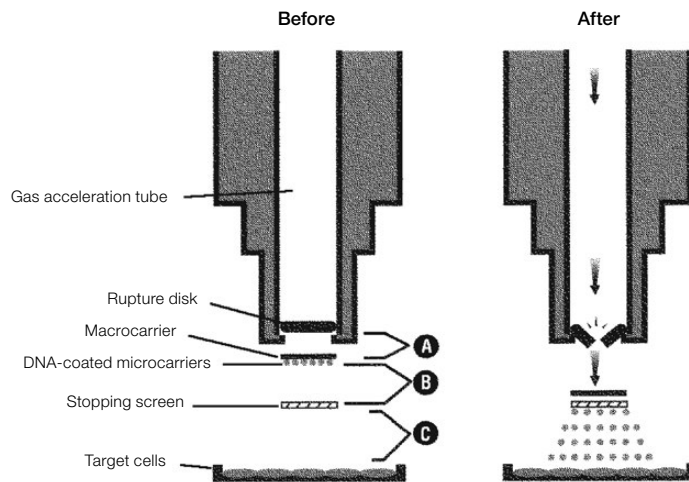


Fig. 1.2. The biolistic bombardment process.

Important Safety Information

Pressurized Helium Safety Information

Caution: Although helium is neither toxic nor flammable, all gases under pressure are potentially dangerous if used improperly. Never use a helium tank with, or attach a tank to, the PDS-1000/He System unless the tank is properly secured. Follow the instructions provided with the helium cylinder from the supplier and those that are applicable for your institution (site safety officer). Bio-Rad has supplied tubing, fittings, a control valve, and a pressure regulator capable of safely handling the high-pressure helium gas used in the biolistic bombardment process. These components have been carefully selected and are the only parts to be used with the PDS-1000/He System.

Power Safety Information

Figure 1.3 shows the serial number certification label, which is found at the rear of the biolistic PDS-1000/He System. This label provides manufacturing data about the instrument and its voltage settings. This instrument and its accessories conform to the IEC and CDRH standards for electrical safety.

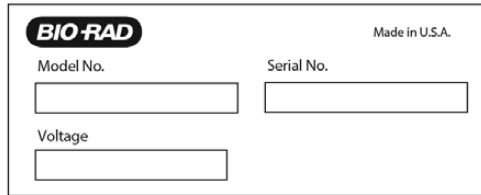


Fig. 1.3. Instrument serial number label on the rear of the instrument.

Requirements for System Operation

Selecting Site for Operation

Prepare a space 61 x 46 x 61 cm (W x L x H) (24 x 18 x 24 in), preferably in a biocontainment hood or other tissue preparation area, near a standard electrical outlet (110 V/60 Hz in the U.S.). Also, allow for placement of the vacuum pump near the site of operation if house vacuum is not used (see Vacuum Supply).

User Supplied Components

Helium Supply

Use only helium gas with the PDS-1000/He System. The low atomic weight of helium permits maximum gas expansion into the bombardment chamber. Thus, sufficient acceleration of the DNA-coated microcarriers is generated to penetrate the target cell membrane.

Obtain a high-pressure (2,400–2,600 psi) tank of high-purity helium to optimize bombardment conditions for the biological system of choice. This allows use of all of the rupture disks (the highest disk has a 1,800 psi rating). Only grade 5 (99.999%) or grade 4.5 (99.995%) helium is to be used, since helium of a lesser grade contains contaminating material, which may obstruct gas flow within the PDS-1000/He System, as well as contaminate the biological sample. Follow all safety instructions provided by the helium supplier for helium tank installation.

The helium pressure regulator (supplied) has a CGA 580, female fitting (standard in the U.S.) for attachment to the user-supplied helium tank. An adaptor to this fitting may be required outside the U.S. Contact your local Bio-Rad office for information on the helium pressure regulator adaptor requirements in your location.

A user-supplied, 1 inch adjustable wrench is required for attachment of the regulator to a helium tank having a capacity of 55 ft³ or greater.

Vacuum Source

The main unit of the PDS-1000/He System must be connected to a vacuum source capable of evacuating the bombardment chamber to a minimum of 5 inches of mercury (5" Hg) for operation. This minimum vacuum requirement is part of the instrument safety system (the chamber door must be sealed for helium pressure to be delivered into the main unit). Tubing is supplied to connect the PDS-1000/He System to vacuum source.

For maximum evacuation capacity, we recommend connecting the PDS-1000/He System to an oil-filled, rotary vane vacuum pump, either single or dual stage, with a pumping speed of 90–150 L/min (3–5 ft³/min). This pumping rate minimizes the time target cells are exposed to vacuum. Oil for the vacuum pump must be supplied by the user. An exhaust mist eliminator on the vacuum pump is also recommended.

The level of the vacuum required in the bombardment chamber depends on the biological system being targeted for transformation. Higher vacuum reduces drag forces on microcarriers during helium-driven acceleration. During the brief bombardment process (less than 1 minute), typical protocols require a vacuum level within the bombardment chamber of 15–29" Hg. Some cells, tissues, and intact plant cells require a high vacuum (up to 28" Hg) for efficient transformation.

House vacuum may be sufficient for bombardment of certain cell types (mammalian cells), but house vacuum pumping rates can fluctuate and vary greatly in overall evacuation capacity (typically 20" Hg, maximum).

Consumables

The 500 Optimization Kit (#1652278) provides the consumables needed for 500 bombardments. It is recommended for users who have yet to determine the optimal conditions for the bombardment of the biological system of interest. The kit contains 0.25 g each of 0.6 μ , 1.0 μ , and 1.6 μ gold microcarriers, 100 each of the seven different rupture disks (ranging from 450 psi to 1,800 psi), 500 macrocarriers, and 500 stopping screens. After optimal conditions are determined, Standard Pressure Kits are available (see Section 7 or the current Bio-Rad catalog for a complete listing).

Additional Laboratory Supplies and Equipment

A vortex mixer is needed for microcarrier preparation. Common laboratory supplies, such as 95% ethanol, pipets, etc. are required, as cited in standard protocols.

Additional Items Available from Bio-Rad

Macrocarrier Holders, set of 5 (#1652322), are included with the biolistic PDS-1000/He System. Additional holders are desirable to facilitate a series of bombardments in one experiment.

Customers outside the continental U.S. and Canada will require a Voltage Converter (#1652259). To use the system, you must locally obtain a cord set that has an IEC 60320/CEE22 connector on one end. This connects to the Voltage Converter and is the type commonly found on computers or televisions. The other end of the power cord will terminate in a plug that will fit the receptacle used in your location. Contact your local Bio-Rad office for more information on this Voltage Converter.

Section 2

Product Description

Packing List

Check the items received with your PDS-1000/He System against the list below. If items are missing, contact your local Bio-Rad office.

- Instruction Manual
- Pressure regulator for helium tank (with in-line 0.45 μm filter)
- 6.0 ft PEEK Plastic Tubing (1/16" OD x .010" ID tubing and fittings), used to connect 3-way helium metering (solenoid) valve to helium pressure regulator

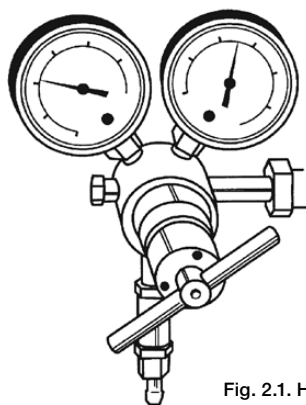


Fig. 2.1. Helium pressure regulator.

- 2.5 ft PEEK Plastic Tubing (1/16" OD x .010" ID tubing and fittings), used to connect 3-way helium metering (solenoid) valve to rear of main unit
- 3-way helium metering (solenoid) valve, with attached cord for power connection to main unit
- Power cord (U.S. and Canada only, 120 V three-prong plug)



Fig. 2.2. 6 ft and 2.5 ft PEEK Tubing.

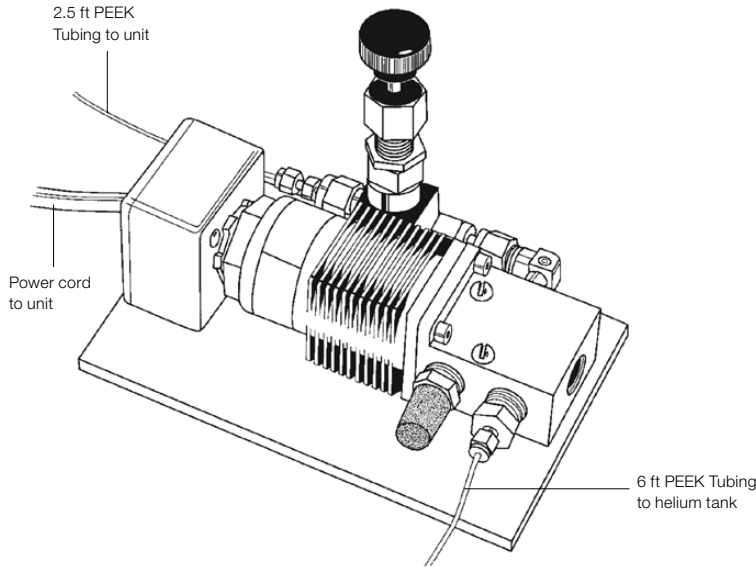


Fig. 2.3. Three-way helium metering (solenoid) valve.

Tools

- 3 hexagonal gap setting tools (1 each 1/8", 1/4", and 3/8"), used to set gap between the bottom of the rupture disk retaining cap and the lid of the microcarrier launch assembly
- One 3/16" hex key (Allen wrench), used for removal of gas acceleration tube (service only)
- One 1/8" hex key, used for releasing set screw in microcarrier launch assembly shelf
- Two 1/4 x 5/16" open-end wrenches, used for connecting small Swagelock Fittings of plastic PEEK Tubing to rear of instrument, external solenoid valve, and helium pressure regulator
- 1 torque wrench for rupture disk retaining cap
- 1 seating tool for macrocarriers



Fig. 2.4. PDS-1000/He tools.

- PDS-1000/He Main Unit (bombardment chamber with control panel and gauges; shipped fully assembled)
- Reinforced PVC vacuum tubing (1/2" ID x 3/4" OD, 5 ft. length and fitting) attached to rear of unit by clamp assembly (centering ring, vacuum hose clamping ring, and nozzle adaptor; see Connecting the PDS-1000/He to a Vacuum Source, in Section 3, for individual components)
- Rupture disk retaining cap (with torque wrench placement holes; Figure 2.6), attached to gas acceleration tube within bombardment chamber

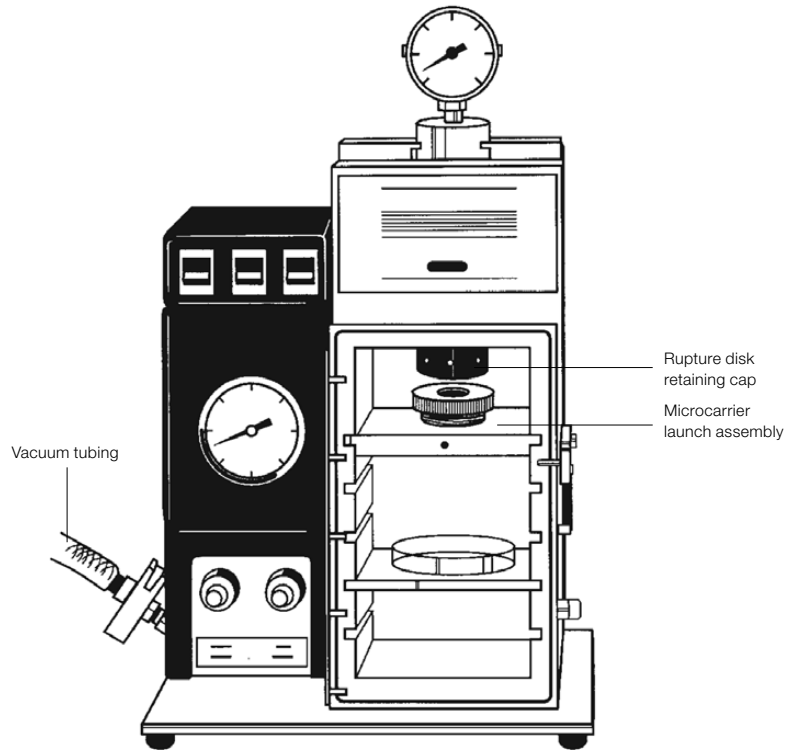


Fig. 2.5. PDS-1000/He Main Unit.

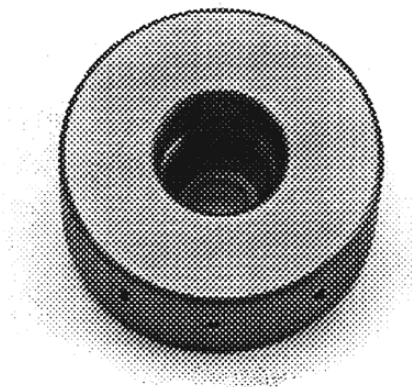


Fig. 2.6. Rupture disk retaining cap.

- Microcarrier launch assembly (shipped fully assembled) consists of the following:
 - Launch assembly shelf with recessed set screw
 - Macrocarrier cover lid
 - Adjustable nest
 - Fixed nest with retaining spring
 - Stopping screen support ring
 - 2 spacer rings, 5 mm height
 - 5 macrocarrier holders, for use within microcarrier launch assembly, after macrocarrier is inserted using macrocarrier insertion tool
- Target plate shelf

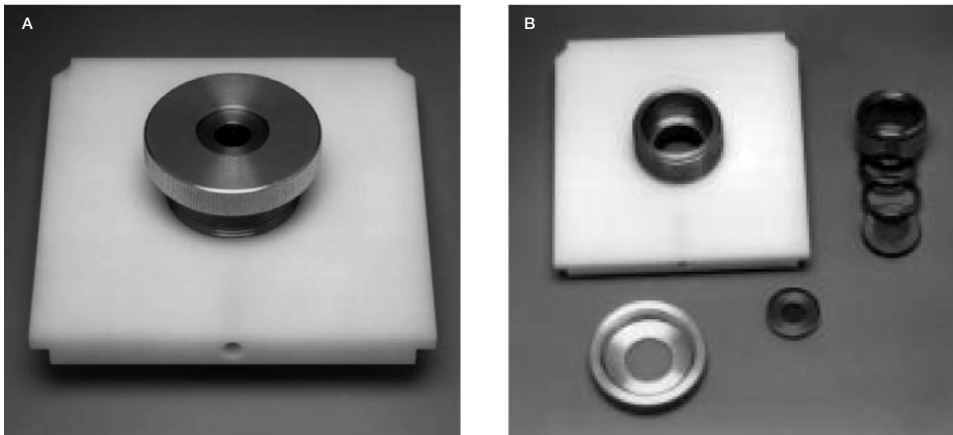


Fig. 2.7. Microcarrier launch assembly (A) and disassembled components (B).

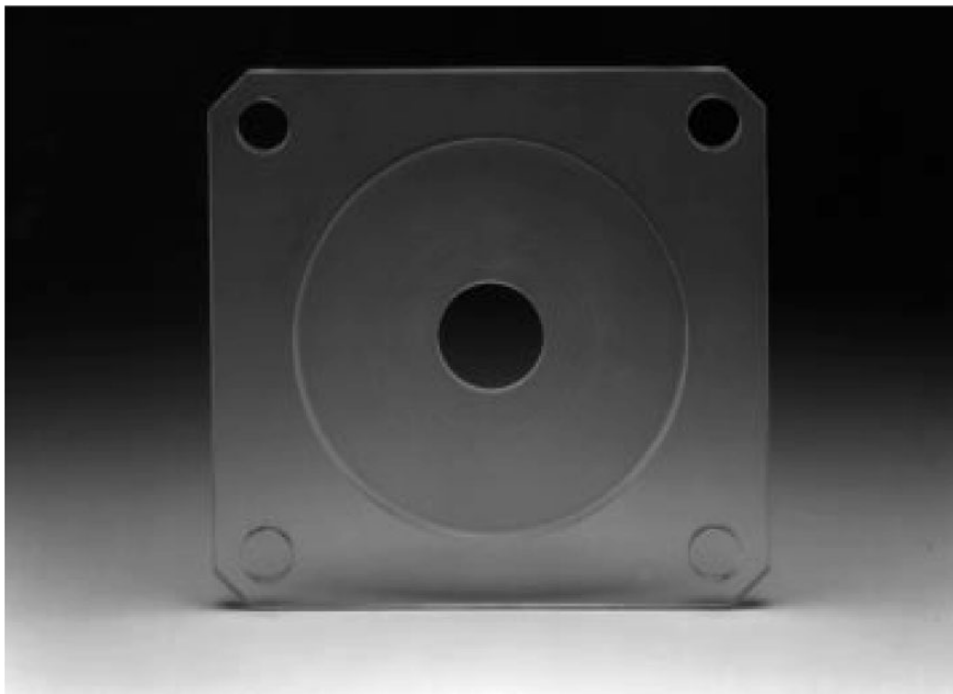


Fig. 2.8. Target plate shelf.

Identification of Unit Controls and Components

The following is a brief description of the operation controls for the PDS-1000/He System (Figure 2.9).

Table 2.1. Unit controls and components.

Controls and Components	Description
Front View, Exterior	
Power switch, ON/OFF	Controls supply of line electrical power to the instrument
Fire switch	Controls flow of helium into gas acceleration tube by activating solenoid valve. Illuminated red when enabled; i.e., when safety interlock is satisfied that at least 5" Hg vacuum is present in chamber Fire switch must be held ON continuously until rupture disk bursts; then release fire switch to stop flow of helium If the fire switch is released before the disk ruptures, the helium is vented via a safety vent in the external 3-way metering (solenoid) valve
Vacuum/vent/hold switch	Controls application of vacuum to bombardment chamber. Vac applies vacuum from line source; Vent releases vacuum using filtered air; Hold maintains vacuum by isolating chamber <i>Bombardments should be performed with this switch in "Hold" position</i>
Vacuum gauge	Indicates level of vacuum in bombardment chamber, in inches of mercury where zero equals ambient atmospheric pressure
Vacuum/vent rate control valves	Regulate rate of application and relief of vacuum in bombardment chamber. Clockwise rotation closes valves
Helium pressure gauge	Indicates helium pressure (in psi) in gas acceleration tube. When solenoid valve is activated by fire switch, the needle in this oil-filled gauge rotates clockwise until rupture disk bursts. <i>Watch this gauge carefully during a bombardment and note the actual rupture pressure</i>
Gas acceleration tube	Helium accumulates within this tube when it is sealed by rupture disk at chamber end of tube. Helium PEEK Tubing connects at top of gas acceleration tube, outside rear of chamber
Bombardment chamber	Holds rupture disk, microcarrier launch assembly, and biological target under vacuum during a bombardment

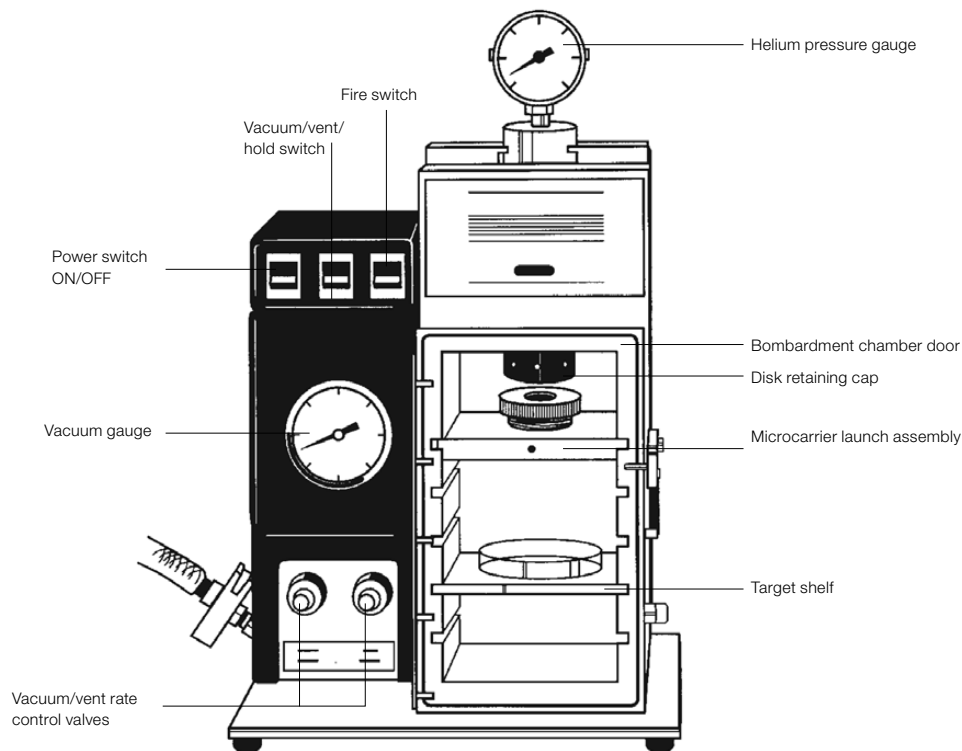


Fig. 2.9. Front view of PDS-1000/He Unit.

Table 2.2. Front view, interior of bombardment chamber.

Bombardment chamber door (with brace)	Controls supply of line electrical power to the instrument
Rupture disk retaining cap	Seals rupture disk against chamber end of gas acceleration tube. This must be tightened securely. The torque wrench is used in the holes in the cap, which are visible in the photo
Microcarrier launch assembly	Holds the DNA/microcarrier preparation on a macrocarrier sheet over the stopping screen in the path of the helium shock wave
Target shelf	Holds the biological target in a petri plate in the path of the accelerated DNA/microcarrier preparation. Particle flight distance is determined by positioning the shelf at one of four levels using slots in the chamber walls

Table 2.3. Rear connections. Refer to Figure 2.10 for a rear perspective view of the unit.

Helium connection to gas tube	Connects top of gas acceleration tube to plastic tubing from solenoid valve, supplying high-pressure helium
Over-pressure relief valve	Opens at 0.5 psi chamber pressure to relieve accumulation of gas. A safety feature. Automatically resets after activation
Vacuum line, chamber to controls	Supplies vacuum to chamber from control valves
Helium metering (solenoid) valve electrical connection	Supplies electric power to 3-way solenoid valve in helium line. Plug the three-pin connector from the solenoid valve into this receptacle
Line cord electrical connection	Supplies electric power to biolistic unit from house line. Plug connector from line cord into this receptacle
Vacuum line, connection to source	Connects unit to house vacuum supply or vacuum pump via fittings included. Feeds into vacuum flow controls

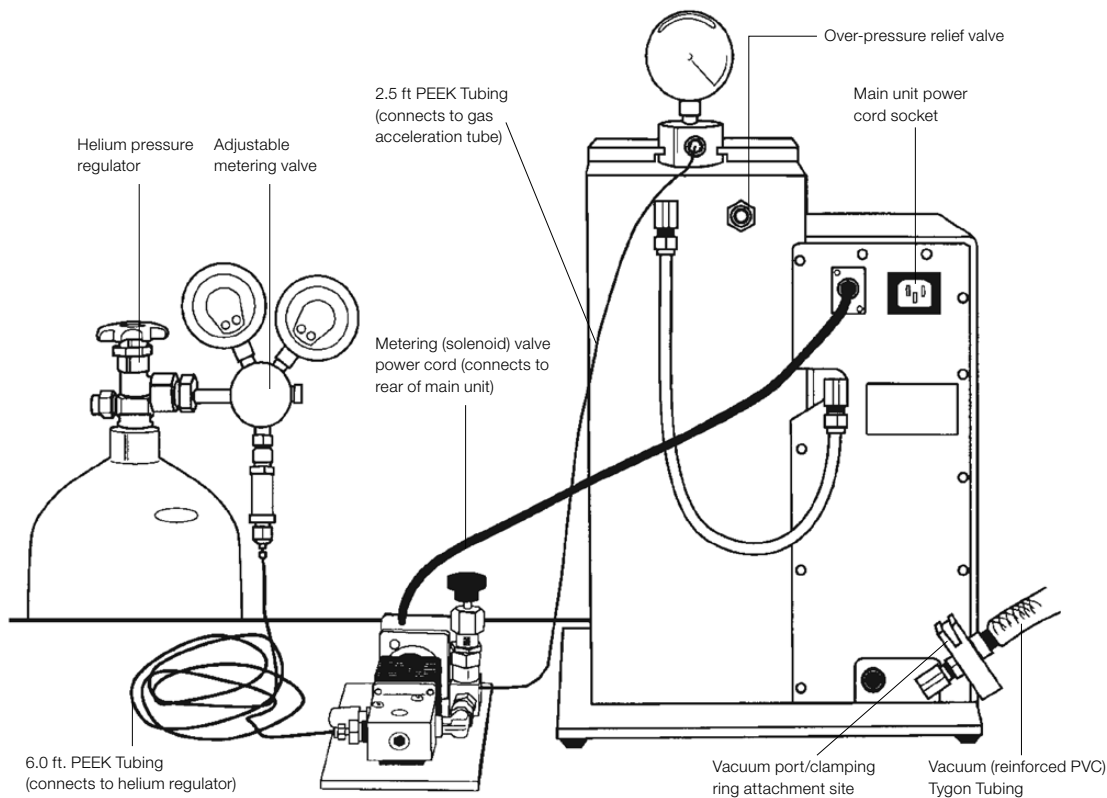


Fig. 2.10. Rear view of main unit, component connection points.

Section 3 Installation

Connecting the PDS-1000/He System to a Helium Source

Refer to Identification of Unit Controls and Components in Section 2 prior to system installation.

Helium Pressure Regulator Installation

Instructions for connecting the helium pressure regulator to a tank of pressurized helium.

Components needed:

- Pressure regulator for helium cylinder (with 0.45 micron in-line filter), provided with unit (Figure 3.1)
- Cylinder of grade 4.5 to 5.0 helium (minimum 99.995% pure); maximum pressure of 2,600 psi, user supplied
- One 1/8" open end wrench or a 10" or 12" adjustable wrench, user supplied

Note: The regulator is intended for use only with helium gas under a maximum of 2,600 psi of pressure. The outlet on pressurized helium cylinders used in the U.S. (maximum 2,600 psi) is compatible with the fitting supplied on the pressure regulator with the PDS-1000/He System (CGA 580, female fitting). Outside of the U.S., contact your local Bio-Rad office for information regarding the proper cylinder/regulator fitting in your area.

1. Secure the cylinder to a wall, post, or other anchored fixture so it will not tip or fall during use.

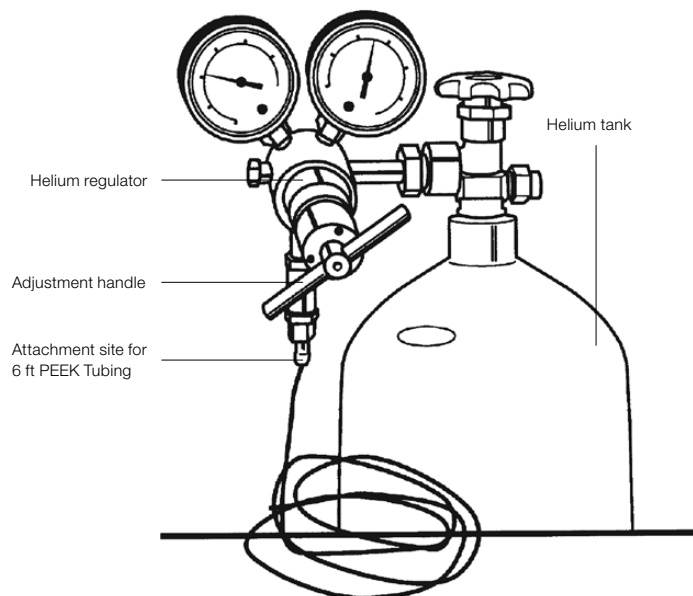


Fig. 3.1. Helium regulator/tank attachment.

2. Inspect the cylinder valve for dirt, dust, oil, grease, or damaged threads. Remove dust and dirt with a clean cloth. Do not attach the regulator if you determine that the valve port is damaged or cannot be cleaned. Inform your gas supplier of this condition and request a replacement cylinder.
3. Clear the valve port of any foreign matter by standing to the side of the cylinder and quickly opening and closing the cylinder valve.
4. Attach the regulator to the cylinder valve and tighten securely with a 1/8" open end wrench or a 10" or 12" adjustable wrench.

Three-Way Metering Valve (Solenoid) Installation

Connecting the helium pressure regulator to the 3-way helium metering (solenoid) valve

Components needed:

- Two 5/16" open end wrenches (Figure 2.4)
- PEEK Tubing, 6 ft (Figure 2.2)
- Solenoid valve (Figure 2.3)

Note: The Swagelock Fittings used on the flexible tubing are easily damaged by overtightening. Turn finger-tight, then use wrench for 1/6 additional turn (60°).

1. Connect one end of the 6 ft tubing assembly to helium pressure regulator (Figure 3.1). Attach the tubing to the in-line 0.45 micron filter at the end of the regulator. Use two 5/16" open end wrenches to tighten the connection.
2. Connect the other end of the 6 ft tubing assembly to the inlet port of the 3-way helium (solenoid) valve assembly. This is the port nearest the metal exhaust filter on the 3-way metering valve. Use two 5/16" open end wrench to tighten the connection (Figures 2.3 and 2.10).

Connecting the 3-way helium metering (solenoid) valve to the PDS-1000/He main unit

Components needed:

- Two 5/16" open end wrenches (Figure 2.4)
- PEEK Tubing, 2.5 ft (Figure 2.2)
- Solenoid valve (Figure 2.3)

Note: The Swagelock Fittings used on the flexible tubing are easily damaged by overtightening. Turn finger-tight, then use the wrench for 1/6 additional turn (60°).

1. Connect one end of the 2.5 ft PEEK Tubing assembly to the outlet port of the 3-way helium metering (solenoid) valve assembly (Figures 2.3 and 2.10). This is the port nearest the adjustable metering valve (black knob). Use two 5/16" open end wrenches to tighten the connection. Do not alter the setting of the metering valve at this time.
2. Connect one end of the 2.5 ft long PEEK Tubing assembly to the PDS-1000/He main unit. This port is located on the gas acceleration tube, below the helium pressure gauge (top, rear of unit). Use a 5/16" open end wrench to tighten the connection (Figure 2.10).

3. Plug the power cord of the 3-way helium metering (solenoid) valve into the rear of the PDS-1000/He main unit (Figure 2.10).

Connecting the PDS-1000/He System to a Vacuum Source

The main unit is shipped with the vacuum hose assembly attached to the port on the rear of the unit via the clamping assembly. Only the connection of the free end of the tubing to a vacuum source is required for connecting the vacuum source to the system.

Choice of Vacuum Source

A vacuum system with a 3 cubic feet per minute (cfm) or 100 L/min pumping capacity is recommended for use with the PDS-1000/He System. This minimizes the total time the target cells/tissue are exposed to reduced atmospheric pressure when inside the bombardment chamber. It is also recommended that the pump have an exhaust mist eliminator.

House or shared source vacuum systems commonly have variable evacuation capacities, from 20 to 23" Hg. When multiple users access a common vacuum system, the actual vacuum pressure will vary, often decreasing to less than 20" Hg per outlet. This is an inadequate vacuum level for yeast and plant cell/tissue bombardment.

Most rotary vacuum pumps will achieve up to 26–28" Hg vacuum. However, the length of time required to attain the desired vacuum depends on the pumping capacity of the vacuum source. The higher the pumping capacity, the less time is required to achieve the desired vacuum. Variation in evacuation time from the bombardment chamber has an unknown effect on target cells and tissue transformation efficiency.

Table 3.1. Pumping rates of various vacuum systems with the PDS-1000/He System.

Vacuum Source Pumping Capacity	Vacuum Level, " Hg	Evacuation Time,* sec
House/common source	20	10–15
	29	Typically not attainable
Pump — 1 cfm or 28 L/min	20	12
	29	45
Pump — 3 cfm or 100 L/min	20	7
	29	25

* Vacuum flow rate valve fully open.

Replacement and Reassembly of Connecting Components

See Figure 3.2 for a diagram outlining component assembly. The clamp consists of an O-ring and a vacuum hose centering ring with male nozzle fitting screwed to it.

1. Place the clamping assembly at one end of the vacuum tubing. The vacuum tubing is PVC-reinforced Tygon Tubing (1/2" ID x 3/4" OD; 5 ft length).
2. The vacuum tubing is shipped with the clamping ring assembly (Figure 3.2) attached to the PDS-1000/He System at the brass vacuum port located on the rear of the unit.
3. Attach the free end of the vacuum hose assembly to the vacuum source.

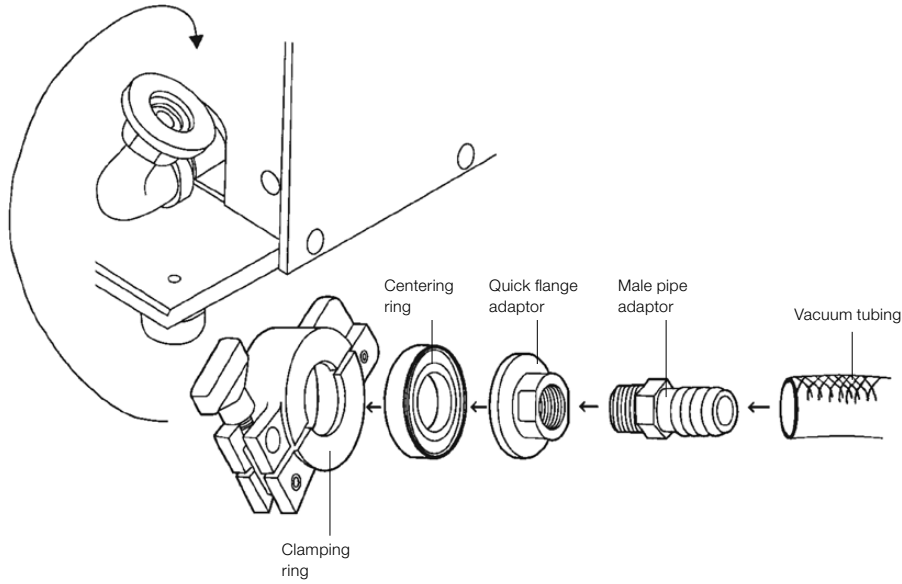


Fig. 3.2. Components used to connect vacuum source to main unit.

Power Cord/Voltage Regulator

The power cord must be plugged directly into a 100–120 V outlet. Plug other end into unit. For use with a 220 or 240 line voltage (users outside the U.S. or Canada), connect a voltage converter to the power cord prior to use (#1652259, see Figure 3.3). You must supply an IEC 60320/CEE22 cord set with a connector on one end that attaches to the voltage converter and a plug on the other end that is compatible with the electrical outlets in your area.



Fig. 3.3. Bio-Rad Voltage Converter.

Section 4

Operation of the PDS-1000/He Instrument

Preparation of System Components Prior to Bombardment

Instrument Preparation

1. Verify that helium tank has 200 psi in excess of desired rupture disk pressure for bombardment.
2. Set gap distance between rupture disk retaining cap and microcarrier launch assembly.

When the rupture disk retaining cap is in place, insert the fully assembled microcarrier launch assembly (with cover lid) inside the bombardment chamber on the highest possible bombardment chamber wall slot (Figure 4.1). The set screw in the white plastic shelf should face outward. Release the set screw on the front of the microcarrier launch assembly with the smaller of the two hex keys provided (Figure 2.4).

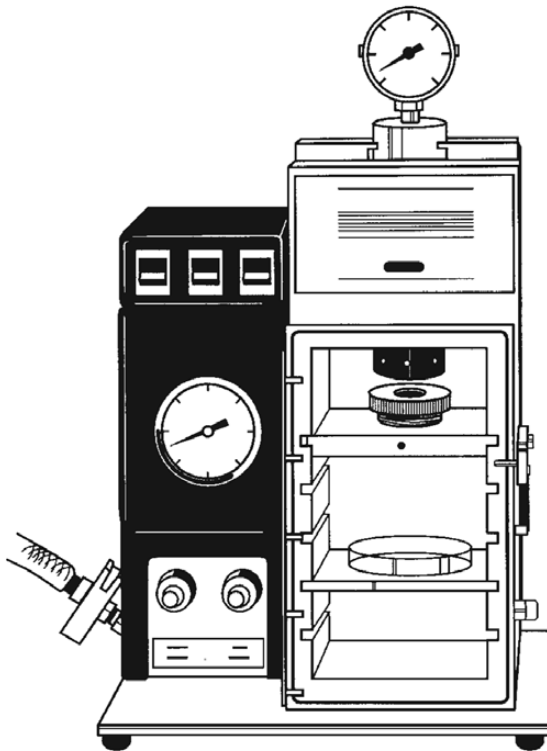


Fig. 4.1. Microcarrier launch assembly and target shelf inside bombardment chamber.

Three hexagonal gap adjustment tools of 1/8, 1/4, and 3/8" have been provided to reproducibly set the gap distance (Figure 2.4). A 1/4" distance between the rupture disk retaining cap and the macrocarrier cover lid is recommended for initial optimization bombardments.

While holding the hexagonal gap adjustment tool against the bottom of the rupture disk retaining cap, turn the adjustable nest until the macrocarrier cover lid touches the gap adjustment tool. The position of the adjustable nest is fixed by tightening the set screw in the white plastic shelf with the hex key.

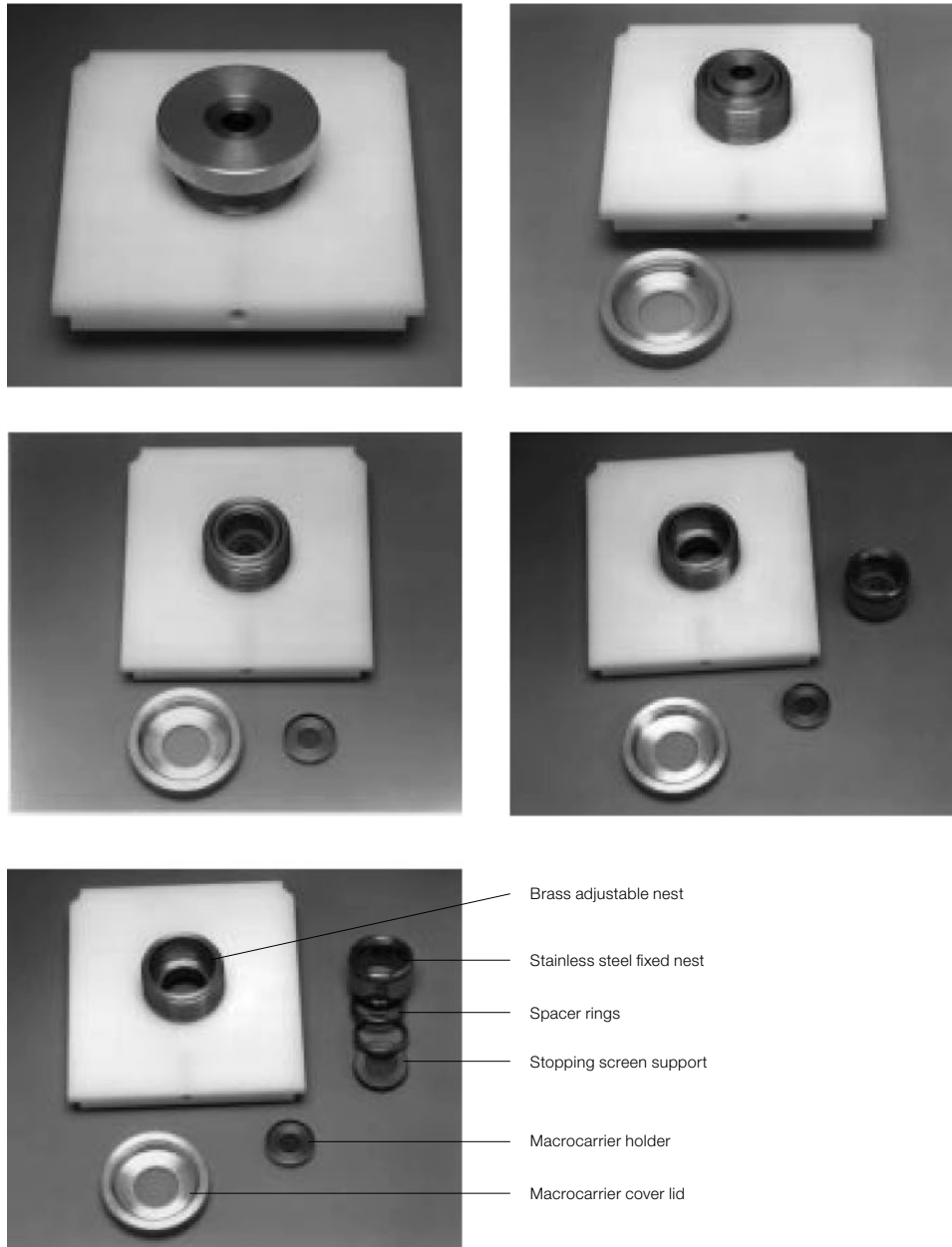


Fig. 4.2. Microcarrier launch assembly adjustment.

3. Prepare the rupture disk retaining cap. After setting the gap between cap and microcarrier launch assembly, wrap the rupture disk retaining cap in aluminum foil and sterilize by autoclaving.
4. Prepare the microcarrier launch assembly. The effect of the gas shock wave on the microcarrier velocities is determined in part by the gap between the rupture disk and the macrocarrier.

The macrocarrier flight distance can be adjusted by varying the position of the stopping screen support and the spacer rings inside the fixed nest (Figure 4.2). The stopping screen support is placed in the middle position at the factory. This is the recommended position when initially optimizing bombardment parameters.

To make this adjustment, remove the microcarrier launch assembly from the PDS-1000/He main unit. Unscrew and remove the macrocarrier cover lid. Disassemble the components of the microcarrier launch assembly by placing the macrocarrier insertion tool into the bottom of the assembly and pushing up. This releases the stainless steel fixed nest from the brass adjustable nest. The two spacer rings and the stopping screen support will fall out from within the fixed nest.

To change the factory-set position to the position for the minimum macrocarrier travel distance (6 mm), invert the fixed nest and insert the stopping screen support inside the fixed nest so that the conical side of the stopping screen support faces down in the final orientation; then insert the two spacer rings (each 5 mm thick). With the fixed nest still inverted, place the microcarrier launch assembly over the fixed nest and seat the fixed nest within the brass adjustable nest.

If greater macrocarrier travel is desired, rearrange the spacer rings and stopping screen support accordingly. The macrocarrier travel distance can be increased in two 5 mm steps, to a maximum of 16 mm.

Sterilize the microcarrier launch assembly by wrapping in aluminum foil and autoclaving. Alternatively, this assembly can be sterilized by wiping with 70% ethanol, followed by drying in a sterile environment.

5. Sterilize the target shelf by wiping with 70% ethanol followed by drying in a sterile environment just prior to use. This part may not be autoclaved.

Consumable Preparation

Several consumables are available for use with the PDS-1000/He System (Figure 4.3).

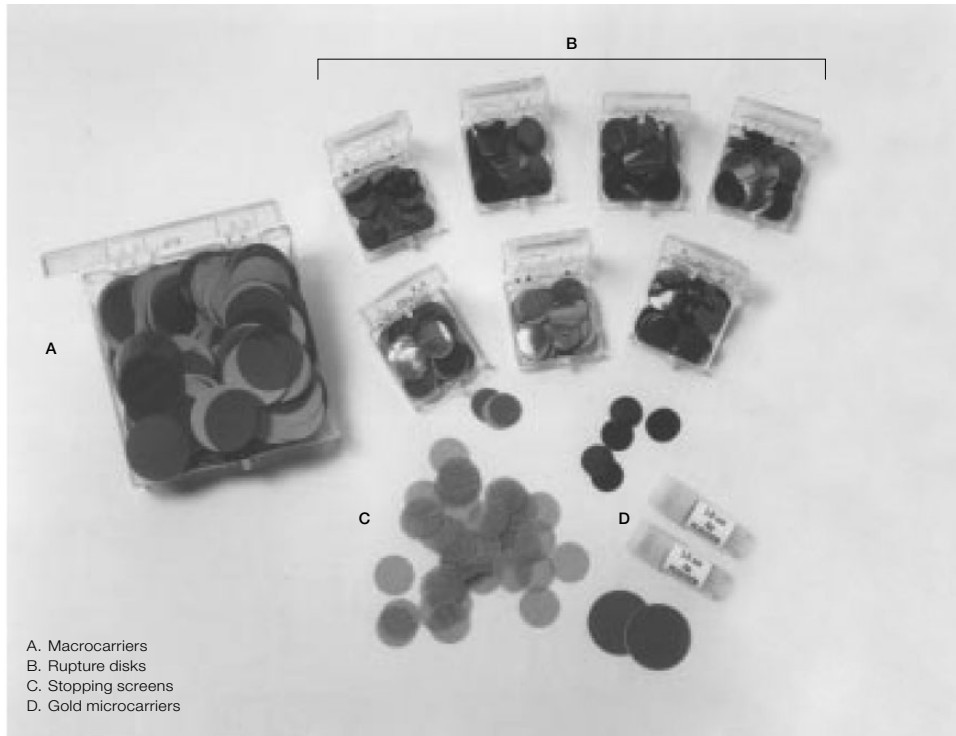


Fig. 4.3. Consumables for the PDS-1000/He instrument. The 0.6 μm gold microcarrier is not pictured.

1. Macrocarriers

Preassemble and presterilize the macrocarrier set in a macrocarrier holder prior to performing sample cell/tissue bombardments.

Macrocarriers are shipped in quantities of 500/box with paper linings between disks. Transfer selected macrocarriers to individual petri dishes for easier handling. Remove the paper lining from between the macrocarriers. Place a macrocarrier into the macrocarrier holder using the seating tool (Figure 4.4). The edge of the macrocarrier should be securely inserted under the lip of the macrocarrier holder. Macrocarrier holders, with macrocarriers already in place, should be sterilized by autoclaving.

2. Rupture disks

Rupture disks (Figure 4.3) are packaged and shipped in quantities of 100/box. Transfer selected rupture disks to individual petri dishes for easier handling. Sterilize rupture disks by briefly dipping them in 70% isopropanol just prior to insertion in the retaining cap. Do not soak for more than a few seconds. Extensive soaking may delaminate the disks, resulting in premature rupture. All disks, with the exception of those rated at 450, 650, and 1,100 psi, are laminated. Autoclaving is not recommended because of potential delamination.

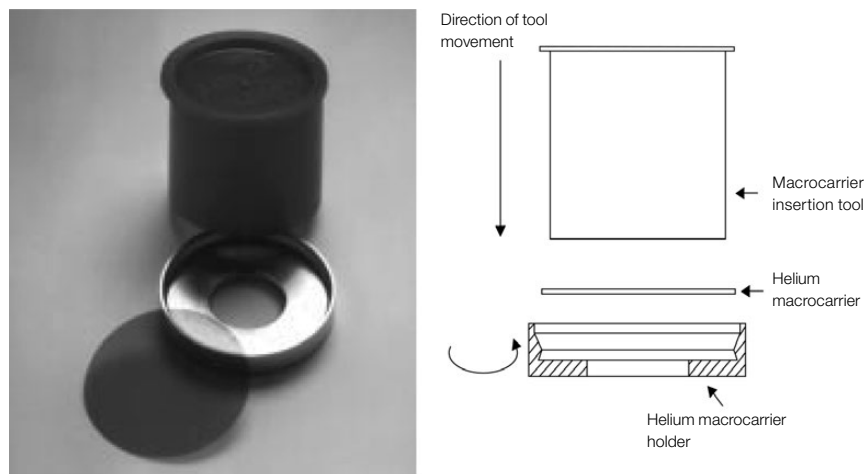


Fig. 4.4. Insertion of macrocarrier into macrocarrier holder with plastic insertion tool.

3. Stopping screens

Transfer selected stopping screens (Figure 4.3) to individual petri dishes for easier handling. Sterilization by autoclaving is recommended. Alternatively, these parts can be sterilized by soaking in 70% ethanol followed by drying in a sterile environment.

4. Microcarriers

The following procedure prepares gold microcarriers for 120 bombardments using 500 μg of the microcarrier per bombardment, based on the method of Sanford, et al. (Sanford et al 1993.)

- a. Weigh out 30 mg of microparticles into a 1.5 ml microfuge tube.
- b. Add 1 ml of 70% ethanol (v/v).
- c. Vortex vigorously for 3–5 min (a platform vortexer is useful).
- d. Allow the particles to soak in 70% ethanol for 15 min.
- e. Pellet the microparticles by spinning for 5 sec in a microcentrifuge.
- f. Remove and discard the supernatant.

Repeat the following wash steps three times:

- Add 1 ml of sterile water
- Vortex vigorously for 1 min
- Allow the particles to settle for 1 min
- Pellet the microparticles by briefly spinning in a microcentrifuge
- Remove and discard the liquid

After the third wash, add 500 μl sterile 50% glycerol to bring the microparticle concentration to 60 mg/ml (assume no loss during preparation).

The microparticles can be stored at room temperature for up to two weeks. Gold aliquots can be stored at 4°C or room temperature.

Store dry gold microcarriers in a dry, nonoxidizing environment to minimize agglomeration.

Coating Washed Microcarriers with DNA

The following procedure is sufficient for six bombardments; if fewer bombardments are needed, adjust the quantities accordingly.

Vortex the microcarriers prepared in 50% glycerol (30 mg/ml) for 5 min on a platform vortexer to resuspend and disrupt agglomerated particles.

When removing aliquots of microcarriers, it is important to continuously vortex the tube containing the microcarriers to maximize uniform sampling. When pipetting aliquots, hold the microcentrifuge tube firmly at the top while continually vortexing the base of the tube.

Remove 50 μ l (3 mg) of microcarriers to a 1.5 ml microcentrifuge tube.

Continuous agitation of the microcarriers is needed for uniform DNA precipitation onto microcarriers. A platform attachment that holds microcentrifuge tubes may be used on your vortex mixer.

While vortexing vigorously, add in order:

- 5 μ l DNA (1 μ g/ μ l)
- 50 μ l 2.5 M CaCl₂
- 20 μ l 0.1 M spermidine (free base, tissue culture grade).

Continue vortexing for 2–3 min.

Allow the microcarriers to settle for 1 min.

Pellet microcarriers by spinning for 2 sec in a microcentrifuge. Remove the liquid and discard.

Add 140 μ l of 70% ethanol (HPLC or spectrophotometric grade). Remove the liquid and discard.

Add 140 μ l of 100% ethanol. Remove the liquid and discard. Add 48 μ l of 100% ethanol.

Gently tap the side of the tube several times and vortex at low speed for 2–3 sec to resuspend the pellet.

Performing a Bombardment

Quick Guide

This summary is repeated in Appendices, Performing a Bombardment, as a tear-out sheet.

Before the Bombardment

1. Select/adjust bombardment parameters for gap distance between rupture disk retaining cap and microcarrier launch assembly. Place stopping screen support in proper position inside fixed nest of microcarrier launch assembly
2. Check helium supply (200 psi over the rated rupture pressure).
3. Clean/sterilize:
 - **Equipment** — rupture disk retaining cap, microcarrier launch assembly
 - **Consumables** — macrocarriers/macrocarrier holders
4. Wash microcarriers and resuspend in 50% glycerol.

5. Coat microcarriers with DNA and load onto sterile macrocarrier/macrocarrier holder on the day of experiment.

Firing the Device

1. Plug power cord from main unit into electrical outlet.
2. Power ON.
3. Sterilize chamber walls with 70% ethanol.
4. Load sterile rupture disk into sterile retaining cap.
5. Secure retaining cap to end of gas acceleration tube (inside, top of bombardment chamber) and tighten with torque wrench.
6. Load macrocarrier and stopping screen into microcarrier launch assembly.
7. Place microcarrier launch assembly and target cells in chamber and close door.
8. Evacuate chamber, hold vacuum at desired level (minimum 5" Hg).
9. Bombard sample by holding Fire button depressed until rupture disk bursts and helium pressure gauge drops to zero.
10. Release Fire button.

After the Bombardment

1. Release vacuum from chamber.
2. Remove target cells from chamber.
3. Unload macrocarrier and stopping screen from microcarrier launch assembly.
4. Unload spent rupture disk.
5. Remove helium pressure from the system (after all experiments for the day have been completed).

Detailed Operation Procedure

Note: We recommend that the first bombardment each day be a “dry run,” with no target cells or microcarriers, to ensure that the system is set up properly and the gas pathway is filled with helium, not air.

1. Power on

With the unit plugged in to the appropriate electrical outlet or voltage converter, turn on the unit by pressing the ON switch. This is the left-most red control panel switch (Figure 2.9).

2. Helium pressure

Confirm that the helium tank pressure regulator is set to 200 psi over the selected rupture disk burst pressure. For example, set the regulator to 1,100 psi when working with a 900 psi rupture disk.

3. Pressurizing the system with helium

Make certain that the helium pressure regulator is properly installed on the helium tank (see Connecting the PDS-1000/He System to a Helium Source, in Section 3).

Close the helium pressure regulator by turning the regulator adjustment screw counterclockwise until the adjusting spring pressure is released.

Release helium into the pressure regulator by carefully and slowly opening the cylinder valve on the helium tank. The cylinder pressure in the tank is indicated on the high-pressure gauge (the gauge closest to cylinder).

Set the desired helium delivery pressure for the rupture disk you are using (measured on gauge on the outside, farthest from cylinder) by turning the regulator adjustment handle clockwise. The pressure should be set to 200 psi above the rupture disk rating.

4. Coating microcarriers with DNA

The day of the scheduled bombardment, coat the microcarriers with DNA. To obtain the best results, use the DNA-coated microcarriers as soon as possible.

5. Loading DNA-coated microcarriers onto a macrocarrier/macrocarrier holder

Each macrocarrier is placed inside a macrocarrier holder and sterilized, as described above. Prior to the application of the DNA-coated microcarriers onto a macrocarrier, prepare a small desiccating chamber for each macrocarrier/macrocarrier holder and place away from excessive vibration.

A small desiccating chamber consists of a sterile 60 mm tissue culture petri dish (with lid) containing CaCl_2 as desiccant in the base of the dish (Figures 4.5 and 4.6). The desiccant is covered with a small piece of filter paper to provide a stable platform for the macrocarrier/macrocarrier holder. The sterile macrocarrier/macrocarrier holder is placed atop the filter paper, with the macrocarrier facing up and the stainless steel holder touching the filter paper.

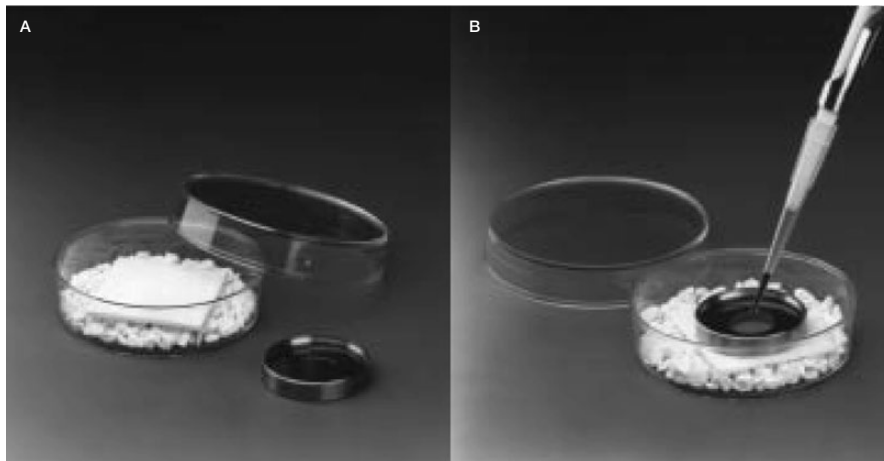


Fig. 4.5A and 4.5B. Loading DNA-coated microcarriers onto a macrocarrier/macrocarrier holder, positioned in small desiccating chamber.

This environment permits rapid access to each macrocarrier and provides a low-humidity environment for the ethanol to uniformly evaporate from the microcarriers. This low humidity, along with a minimum of vibration during evaporation, minimizes microcarrier agglomeration.

For each macrocarrier, remove 6 μl aliquots (approximately 500 μg) of microcarriers and spread evenly over the central 1 cm of the macrocarrier using a pipet tip. Pipet from a continuously vortexed tube and rapidly apply suspended microcarriers to the macrocarrier, as microcarriers quickly settle out from the ethanol solution in the tube or even in the pipet tip.

Cover the culture dish immediately after application of the microcarrier suspension to the macrocarrier. The ethanol should evaporate within 10 min to leave the DNA-coated microcarriers adhering to the macrocarrier. The loaded macrocarriers should be used within 2 hr.

6. Cleaning chamber walls

Clean the PDS-1000/He chamber as desired with 70% ethanol. Allow time for drying. Do not autoclave or flame sterilize the PDS-1000/He System.

7. Loading the rupture disk

Unscrew the rupture disk retaining cap from the gas acceleration tube from within the bombardment chamber (Figure 4.6) or unwrap cap from sterile wrapping.

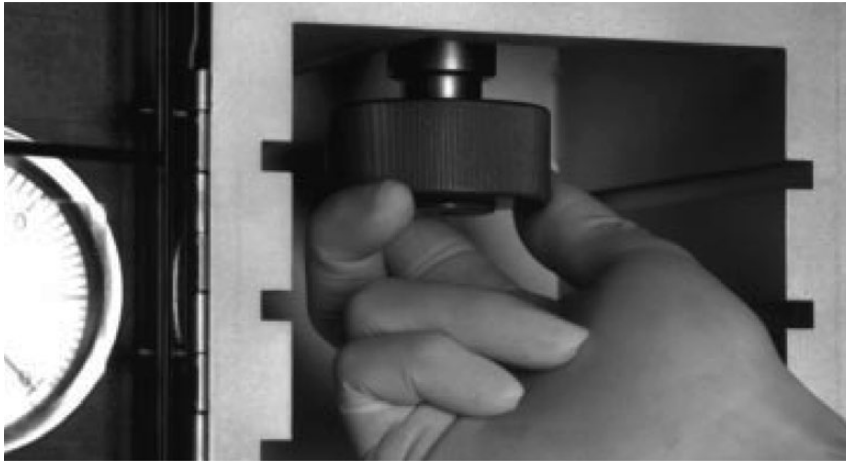


Fig. 4.6. Removing/mounting of rupture disk retaining cap onto end of gas acceleration tube inside bombardment chamber.

Select rupture disk of desired burst pressure. Handle all rupture disks with sterile forceps. Grease, fingerprints, or even powder from plastic gloves on the rupture disk may prevent a tight seal from forming within the retaining cap.



Fig. 4.7. Rupture disk insertion into recess of retaining cap.

Immediately before placing a rupture disk in the retaining cap, briefly wet the rupture disk in isopropanol (do not soak the disk for an extended period of time or the disk may delaminate). Loading the rupture disk while wet forms a liquid gasket when the cap is screwed onto the gas acceleration tube, and thereby reduces the failure rate of the rupture disk. Place the rupture disk in the recess of the rupture disk retaining cap (Figure 4.7).

Screw the rupture disk retaining cap onto the gas acceleration tube using a left-to-right motion. Never tighten the rupture disk retaining cap without a rupture disk in place or scratching and deformation of the two metal surfaces will occur and cause helium to leak when a rupture disk is pressurized.

The retaining cap is tightened to a torque of 60 in/lb with the retaining cap torque wrench. To use the torque wrench, insert the short end of the metal rod into an accessible hole in the retaining cap. Push the long end of the rod to the right only until it touches the inner surface of the black tube (Figure 4.8). If the retaining cap is not tightened sufficiently, the rupture disk may slip out of place as the gas acceleration tube fills with helium. Test fire once to fill gas tubing with helium.



Fig. 4.8. Proper torque applied to retaining cap with torque wrench.

8. Microcarrier launch assembly

Unscrew the macrocarrier cover lid from the assembly. Place a sterile stopping screen on the stopping screen support (Figure 4.9). Note: Never operate the PDS-1000/He Instrument without a stopping screen in place. The target sample will be destroyed from the uninterrupted acceleration of the macrocarrier by the helium shock wave.

Install the macrocarrier/macrocarrier holder on the top rim of the fixed nest (Figure 4.10). The dried microcarriers should be facing down, toward the stopping screen. Replace the macrocarrier cover lid on the assembly and turn clockwise until snug. Do not overtighten.

Place the microcarrier launch assembly in the top slot inside the bombardment chamber (Figures 4.11 and 5.1).

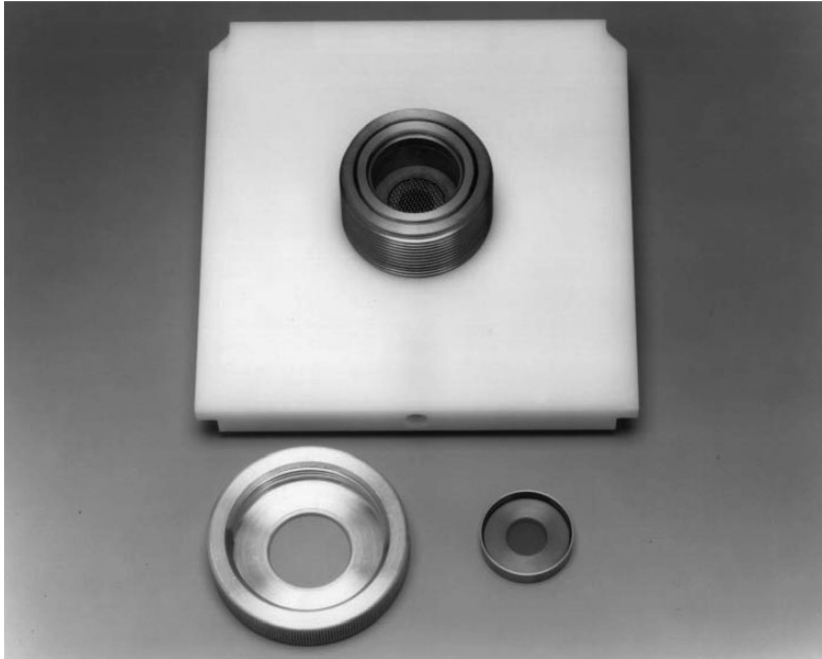


Fig. 4.9. Placement of stopping screen inside fixed nest with macrocarrier and cover lid removed.



Fig. 4.10. Removal/replacement of macrocarrier cover lid with assembled fixed nest. Macrocarrier holder (with macrocarrier properly inserted) is inverted and placed atop fixed nest.

9. Target cells/tissue placement in chamber

Place the target shelf at the desired level inside the bombardment chamber. Place the sample (usually contained within a petri dish) on the target shelf. Close and latch the sample chamber door.

10. Chamber evacuation/hold

Turn on the vacuum source. Set the vacuum switch on the PDS-1000/He System (middle red control switch, Figure 4.11) to the **VAC** position. Evacuate the sample chamber to the desired level, at least 5" Hg. The right-most red control switch (the **FIRE** switch) will be illuminated when the minimum vacuum is achieved.

When the desired vacuum level is reached, hold the chamber vacuum at that level by quickly pressing the vacuum control switch through the middle **VENT** position to the bottom **HOLD** position.

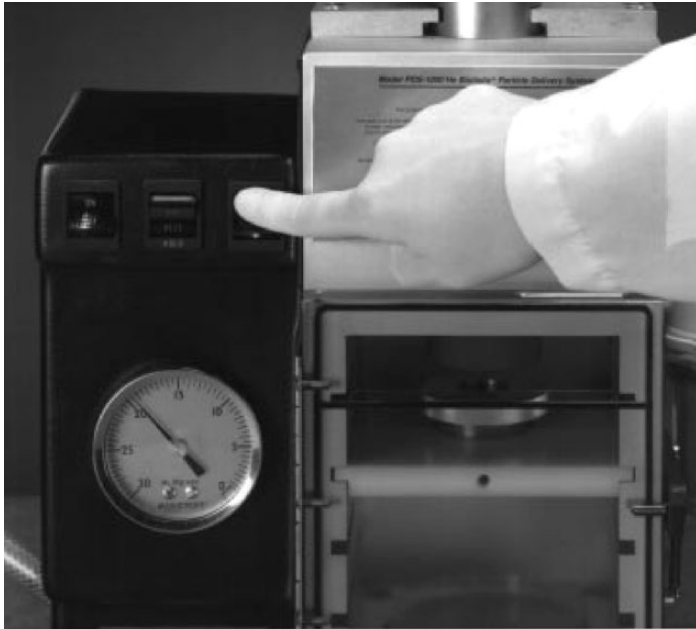


Fig. 4.11. Sample bombardment: power switch ON, vacuum switch on HOLD position, and user continuously pressing FIRE button.

11. Bombard the sample

With the vacuum level in the bombardment chamber stabilized, press and hold the FIRE switch to allow helium pressure to build inside the gas acceleration tube that is sealed by a selected rupture disk (Figure 4.11).

Estimate rupture disk burst pressure by observing the helium pressure gauge at the top of the acceleration tube. A small pop will be heard when the rupture disk bursts. The rupture disk should burst within 10% of the indicated rupture pressure and within 11–13 sec.

Release the **FIRE** switch immediately after the disk ruptures to avoid wasting helium. Releasing the **FIRE** switch prior to disk rupture will vent the gas acceleration tube via the 3-way helium metering (solenoid) valve.

Note: Variation in the burst pressure indicated on the helium pressure gauge (on the top of the unit) from the rated rupture disk pressure may be observed if the gas acceleration tube fill rate is improperly set. See Solenoid Valve Adjustment, in Section 5.

12. Release vacuum from chamber

Release the vacuum in the sample chamber by setting the **VACUUM** switch to the middle **VENT** position.

13. Target cells removal from chamber

After vacuum is released, the vacuum gauge should read 0" Hg of vacuum. Open the sample chamber door. Remove the sample and treat as appropriate.

14. Macrocarrier and stopping screen removal from microcarrier launch assembly

Remove the microcarrier launch assembly. Unscrew the lid and remove the macrocarrier holder. Discard the used macrocarrier and stopping screen (Figure 4.12).

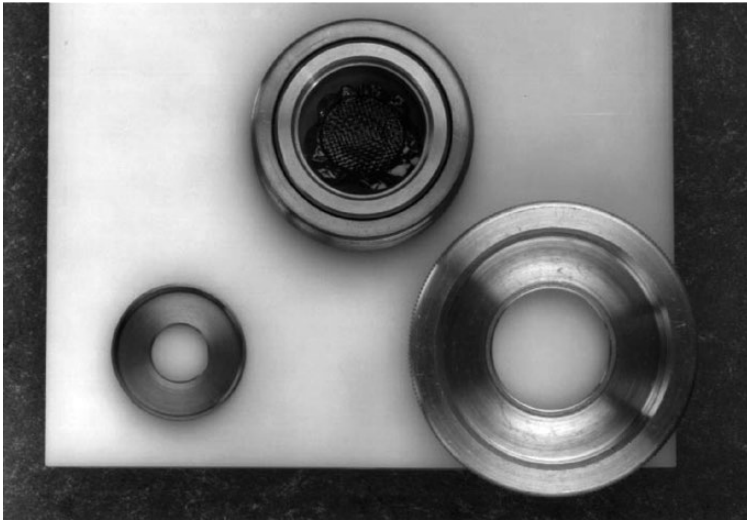


Fig. 4.12. View of used macrocarrier and stopping screen within disassembled microcarrier launch assembly after a bombardment.

15. Removal of spent rupture disk

Unscrew the rupture disk retaining cap from the gas acceleration tube. Remove the remains of the rupture disk (Figure 4.13). The next bombardment may now be performed (from step 7).

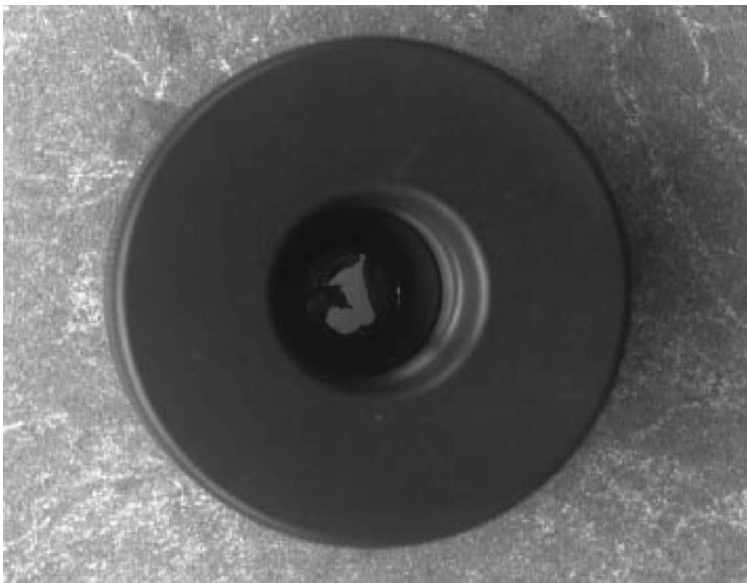


Fig. 4.13. View of spent rupture disk within the retaining cap.

Removal of Residual Helium Pressure – Shutdown

After completing all bombardment(s), remove the helium pressure from the PDS-1000/He System and close the helium cylinder valve. Perform the following steps to remove helium pressure from the system.

1. Close the helium cylinder valve and chamber door.
2. With at least 5" Hg of vacuum in bombardment chamber, remove residual line pressure from the regulator, solenoid, and PEEK Tubing by activating the **FIRE** button.
3. Release the **FIRE** button on the apparatus, and remove all tension on the pressure adjustment screw of the helium regulator by turning counterclockwise until it turns freely.
4. Vent any residual vacuum from the bombardment chamber by setting the vacuum switch to the **VENT** position.

Section 5

Selection and Adjustment of System Bombardment Parameters

Overview – Matrix of Variables

Factors affecting bombardment efficiency are numerous, and they interact in complex ways. All possible variables cannot be addressed in a single large factorial experiment. Instead, most users will find it sufficient to prioritize variables by the magnitude of their effects, optimizing them individually, and then testing their interactions on a more limited scale. Table 5.1 shows some examples of conditions used in published protocols for a variety of target materials.

Table 5.1. Cell types, settings, and conditions.

Cell Type	Growth Phase	Cell Density	Osmoticum	Vacuum, " Hg	Target Distance, cm	Helium Pressure, psi	Particle Size
Bacteria	Late log to early stationary	10 ⁸ –10 ⁹ per 100 mm plate	0.75 M sorbitol	29	6	1,100	M5 tungsten
Yeast	Early stationary	10 ⁸ –10 ⁹ per 100 mm plate	0.75 M sorbitol and 0.75 M manitol	28	6	1,300	0.6 µm gold
Algae	Log	10 ⁸ –10 ⁹ per 100 mm plate	—	29	6	1,300	0.6 µm gold
Plant	—	10 explants per 100 mm plate	None	28	6	1,300	1.0 µm gold
• embryos	—	10 explants per 100 mm plate	None	28	6	1,300	1.0 µm gold
• callus or cell culture	Log	0.75 ml packed cell volume	None	28	9	1,300	1.0 µm gold
Subcellular organelles	Mid-log	5 x 10 ⁷ per 100 mm plate	None	28	6	1,300	0.6 µm gold
Animal	Log	50–80% confluent on 35 mm plates	None	15	3	1,100	1.6 µm gold
• tissue culture	Log	50–80% confluent on 35 mm plates	None	15	3	1,100	1.6 µm gold
• tissue sections	1 hr–4 days post-excision	400 µm sections	None	25	9	1,100	1.6 µm gold

Vacuum Level in Bombardment Chamber

The vacuum in the bombardment chamber reduces the frictional drag of the microcarriers as they are accelerated toward the target cells. The unit should be connected to a vacuum system (see previous section) that can evacuate the bombardment chamber to 28–29" Hg in less than 30 seconds. This level of vacuum is useful for most plant and microbial cells/tissues. Mammalian cells/tissue should be bombarded at approximately 15" Hg.

Helium gas enters the evacuated bombardment chamber once the rupture disk bursts. The use of low molecular weight helium minimizes the deceleration of the microcarriers as they pass through helium and also reduces the force of the gas shock wave that hits the target cells. This reduced impact will help minimize target tissue damage.

Helium Pressure/Rupture Disk Selection

Each of the seven different rupture disks available ruptures at a specific pressure, ranging in rating from 450 to 1,800 psi. The rupture pressure determines the power of the shock wave entering the bombardment chamber. Increasing helium pressure will increase particle acceleration and subsequent target tissue penetration by the DNA-coated microcarriers. Since the shock wave or resulting acoustic wave may cause damage to the target cells or tissue, use the lowest helium pressure that gives high transformation efficiency.

Table 5.2. Suggested starting helium pressure conditions for optimizing various biological systems.

Cell Type	Rupture Disk Pressure, psi
Bacteria	1,100
Fungi	1,300
Yeast	1,300
Plant cells/tissue	1,100
Mammalian cells	1,100

Solenoid Valve Adjustment

A factory preset metering valve (black knob) on the 3-way helium metering (solenoid) valve assembly controls the fill rate of the gas acceleration tube. The proper fill rate is set at the factory for a 1,550 psi rupture disk to burst within 12–15 seconds. A more rapid fill rate may result in what appears to be a lower burst pressure, due to a lag in needle movement in the oil-filled gauge. The metering valve (black knob) may be adjusted if desired. Adjust knob in small increments: a clockwise rotation of the black knob will lengthen the fill rate.

Vacuum Flow Rate Control Valves

Both valves are set at the factory to be fully open, for maximum flow rates. To decrease the rates, turn the valves (knob) clockwise until the desired evacuation and/or vent rate is achieved (see Figure 2.9).

Distance between Rupture Disk and Macrocarrier

The effect of the gas shock wave on microcarrier velocities is determined in part by the gap between the rupture disk and the macrocarrier. The smaller the distance, the more powerful the effect of the gas shock wave on macrocarrier acceleration.

Three hexagonal gap adjustment tools of 1/8, 1/4, and 3/8" have been provided to reproducibly set the gap distance (Figure 2.4). A 1/4" distance between the rupture disk retaining cap and the macrocarrier cover lid is recommended when initially optimizing bombardment parameters.

Distance between Macrocarrier and Stopping Screen

Macrocarrier flight instability increases with greater travel distance. Therefore, the shortest travel distance is recommended. This is achieved by having the stopping screen support placed above both spacer rings inside the fixed nest when initially optimizing bombardment conditions. The macrocarrier flight distance can be adjusted by varying the position of the stopping screen support, as positioned by the spacer rings inside the fixed nest of the microcarrier launch assembly. The travel distance can be increased to a maximum of 16 mm in steps of 5 mm by varying spacer ring placement.

Distance between Stopping Screen and Target Shelf (Microcarrier Flight Distance)

One of the most important parameters to optimize is target shelf placement within the bombardment chamber. This placement directly affects the distance that the microcarriers travel to the target cells for microcarrier penetration and transformation.

Four target shelf levels are available in the bombardment chamber: level 1 = 3 cm, level 2 = 6 cm, level 3 = 9 cm, and level 4 = 12 cm below the stopping screen (Figure 5.1).

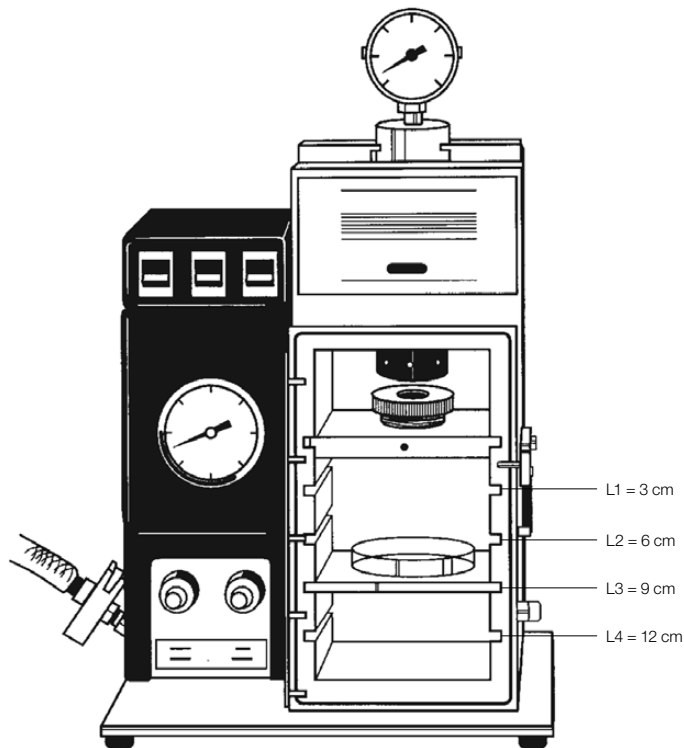


Fig. 5.1. Target shelf placement and corresponding target distances.

Microcarrier Selection

Types of Microcarrier

Two types of microcarriers can be used with the PDS-1000/He System, gold and tungsten. The density of these microcarriers is sufficient to penetrate a wide variety of cell and tissue types using the biolistic PDS-1000/He Acceleration System.

Gold microcarriers are available in three sizes. Gold particles are more spherical and are more uniform in size within a given sample than tungsten microcarriers. Gold is biologically inert, but the DNA coating of this type of microcarrier is subject to more variation during precipitation than tungsten.

Tungsten particles can also be used with the PDS-1000/He System. They are more irregular in shape and size variation within a given particle size and are much wider than gold. The size variation makes optimization using a certain particle size more difficult. Tungsten may be toxic to certain cell types and may oxidize to alter DNA binding.

Microcarrier Size

The velocity of the chosen microcarrier increases with larger particle size. Increased microcarrier velocity has a major effect on particle penetration into target cells/tissue and thus transformation efficiency.

Table 5.3. Recommended starting particle size/type for bombardment of various cell types.

Cell Type	Particle Size
Bacteria	0.7 μm (M5) tungsten
Yeast	0.6 μm gold
Algae	0.6 μm gold
Plant cells/tissue	1.0 μm gold
Animal cell cultures	1.6 μm gold
Subcellular organelles	0.6 μm gold

Table 5.1 and bulletins 1688 and 2015 give a detailed discussion of optimization parameters.

Preparation of Biological Material for Bombardment

Plant Cells

Many factors contribute to optimum performance in any system for plant gene transfer. Transformation procedures employing microcarrier bombardment are no exception, but the improved PDS-1000/He System lends itself to rapid and easy modification to tailor its performance to a particular application. Microcarrier bombardment also allows exploitation of a very broad range of explant tissue types. This permits optimization of a crucial aspect of the transformation system, which is difficult to alter when transforming by other means.

The following is a discussion of some of the critical parameters in the microcarrier transformation system, together with a brief examination of explant characteristics that affect transformation efficiency. This is not intended as a “how to” manual. Because of the extreme diversity of applications possible with the system and the various types of tissues that can be employed, it is not possible to produce a universal protocol for all applications. Rather, this discussion is intended to aid the new user to design experiments.

Choice of Explant Tissue

Unlike most other procedures for DNA transfer, microcarrier bombardment places few constraints on the types of tissues that can be treated. DNA can be delivered into any cell or tissue that can be exposed sufficiently to allow particle penetration. Thus more attention can be given to issues such as the tissue's ability to regenerate plants, or its physiological suitability for gene regulation studies. There are, however, several aspects of the explant that contribute to the efficiency of the system.

An ideal tissue for microcarrier bombardment could perhaps be described as a cell monolayer, all cells of which are capable of expressing introduced genes, and all of which could, independently, divide and differentiate into functional plants. Microcarrier bombardment of such a tissue will likely be efficient because cells can be spread over a large area, thus efficiently capturing a large proportion of the particles delivered in each bombardment. Particles also have a high probability of penetrating such cells because they are not covered by overlying cell layers which would very likely intercept some of the particles delivered by each bombardment. Absolute transient expression levels are likely to be high, because of the number of cells affected. The low profile of cell clusters reduces the probability that the tissue will be moved about by the gas released in the shock wave during treatment. Such a tissue would also lend itself well to selection with antibiotics or herbicides, in part, at least, because it would have a favorable geometry for the establishment of a uniform concentration of the selective agent throughout the culture.

Some kinds of embryogenic suspension cultures approach this ideal. In fact, it is this type of tissue that has allowed great success in the microcarrier-mediated transformation of recalcitrant species such as maize (Gordon-Kamm et al. 1990). While most explants differ from this ideal in substantive ways, it is useful to keep these characteristics in mind when choosing and preparing tissues for bombardment, whether taken from established cultures or recently removed from donor plants. For example, tissue damage is typically most severe within the central area of bombardment where particle density is greatest. Cultured materials can usually be arranged to minimize this effect by placing them in a circle around the central area of bombardment. Even though some of the DNA-bearing particles fail to enter cells, this will often increase the absolute number of transformation events recovered from a single bombardment by reducing tissue damage. Similarly, explants from donor plants should be arrayed so as to capture as large a portion of particles as possible. Newly acquired explants can be dissected so that the cell types of interest are exposed as much as possible. Minimize the height of the tissue above the surface of its support medium to reduce losses during treatment, which result from tissues with high profiles being blown about during bombardment. Finally, it may be useful to reduce the size of each explant or tissue mass to more closely approximate the geometry of suspension culture cells and enhance the uniformity of selection.

Most tissues appear to vary with regard to their ability to express introduced genes following bombardment. This effect is frequently observed as fluctuations in the frequency or intensity of transient expression of β -glucuronidase (GUS) constructions among different tissues, or even among cells in different areas of a single explant. It can be caused by uneven particle distribution during preparation of macrocarriers and also by heterogeneity of the tissue itself. Variation in expression associated with an explant's physiological age is not uncommon in intact explants. Although the causes of these phenomena are unknown, it is important to realize that such effects exist and can be confounding in experiments designed to choose tissues for transformation experiments and to optimize bombardment parameters.

Optimization of Bombardment Parameters — Transient Expression vs. Stable Transformation

The goal of many microcarrier bombardment experiments is to develop protocols for efficient production of transgenic plants. The frequency at which stable transformation events are recovered is the ultimate criterion by which these protocols must be measured. However, measurement of transient expression (that is, expression of newly introduced DNA sequences shortly after bombardment) provides rapid feedback and can be invaluable in determining effective bombardment parameters for specific tissues.

A precise quantitative relationship between the level or frequency of transient expression observed in a bombarded tissue, and the frequency at which stable transformation of cells occurs within that tissue may not exist. Transient expression is, however, a very useful indicator of the efficiency of DNA delivery, and can help to define conditions required to deliver DNA into specific cell layers. In this way, measurement of transient expression allows rapid determination of a set of parameters which permit delivery of DNA into tissues of interest, and provides critical information about which parameters have the greatest effect.

Choice of reporter gene and assay procedures for optimization of bombardment must be determined by the particular question to be addressed. Histochemical assay of β -glucuronidase expression in bombarded cells is ideal for determining which cell types in a heterogeneous tissue are capable of expressing the introduced sequence. This procedure is not useful, however, for determining the absolute expression levels produced by a particular set of bombardment parameters. For such measurements, it is more appropriate to use an extractive assay procedure in which expression level can be measured, for example, as a function of enzyme activity per unit of total protein. Several excellent markers (and associated assay procedures) are available for this type of study. These include luciferase and β -glucuronidase measured by a fluorometric process, and neomycin phosphotransferase, for which a commercial ELISA assay is available.

Yeast/Bacteria/Microbes

Cells that grow in suspension should be grown to late log to early stationary phase in media under the appropriate conditions. Pellet the cells and resuspend in sterile water. Estimate the cell density by determining the absorbance (for yeast, 1 OD₆₀₀ is approximately 2.8×10^9 cells/ml; for *E. coli*, 1 OD₆₀₀ is approximately 1.0×10^9 cells/ml). Spread 1×10^8 yeast or 2×10^9 *E. coli* on 100 mm petri plates containing the appropriate media. In addition to the necessary supplements, the media for growing yeast should contain 0.75 M mannitol and 0.75 M sorbitol, while media for growing bacteria should contain 0.75 M sorbitol. The plates may be allowed to air dry briefly, but should be used within 1 or 2 hr. Transformation efficiency decreases if the plates are allowed to sit even for several hours.

Mammalian Cells

Cells in Culture

Both primary and established animal cell lines are transformable using particle bombardment. Because optimum bombardment conditions are usually found when mammalian cells are placed at the top level of the bombardment chamber, the cells can be inoculated into 35 mm tissue culture plates. Only if cells are to be bombarded at Target Level 2 or farther should larger plates be used.

Cells that grow as attached cultures should be inoculated into tissue culture plates one day prior to bombardment so that they are in log phase and 50–80% confluent at the time of transformation. Cells that grow as suspension cultures should be inoculated from log phase cultures onto plates that have been aseptically coated by adsorption with Cell-Tak (Thermo Fisher Scientific, Waltham, MA) according to the manufacturer's instructions. Allow cells to attach for 1–2 hr. They should be approximately 75% confluent at the time of bombardment.

Aspirate tissue culture media immediately prior to bombardment so that the cells are directly exposed to air. Place cells to be bombarded in the bombardment chamber, centered under the stopping screen, and close the chamber door. Evacuate the chamber to the proper vacuum and bombard the cells. After bombardment, add fresh tissue culture media to the cells and incubate for the appropriate time.

Organ Cultures

Animal organs may be transformed by particle bombardment 1 hour to several days after preparation. Tissues should be in sections approximately 400 μm thick and no more than 1 cm square. The tissue may be grown in cell culture inserts containing a permeable membrane. The tissue section (in the cell culture insert) is placed in a petri plate containing 1% agar and then transferred to the bombardment chamber. After bombardment, the insert containing the tissue can be transferred back to the growth chamber.

Section 6 Troubleshooting

Rupture Disk Bursts at Incorrect Pressure

Probable Cause	Action
1. The helium flow rate is too fast	Note the helium burst pressure by observing the helium pressure gauge at the top of the main unit. A factory preset metering valve (black knob) on the external 3-way helium metering (solenoid) valve assembly controls the fill rate of the gas acceleration tube. The proper fill rate is 12–15 seconds to burst a rupture disk when the helium pressure is set 200 psi over the rated rupture pressure. A faster fill rate may result in what appears to be a lower burst pressure due to a lag in needle movement in the oil-filled gauge. The metering valve (black knob) may be adjusted. Adjust knob in small increments: a clockwise rotation of the black knob will lengthen the fill rate
2. An improper disk was inserted or two disks are sandwiched together	Replace with a single rupture disk of the correct psi
3. The retaining cap was not securely tightened	This is frequently the reason when a disk slips out of the retaining cap. It is important to screw the cap onto the end of the gas acceleration tube by hand and then use the torque wrench
4. The rupture disks are not clean	Use tweezers to handle the disks after briefly rinsing in isopropanol. Do not handle with bare or gloved hands (grease and powder source)
5. The sealing edge (inside bombardment chamber) of the gas acceleration tube is not clean	Occasionally adhesive from the (laminated) rupture disks may build up on the edge of this part. The buildup can be cleaned by using a mild, nonabrasive detergent, like dish soap
6. The sealing edge of the gas acceleration tube is damaged	The edge of this part is susceptible to physical damage if the retaining cap is tightened without a rupture disk in place. If damage has occurred such that the disk will not rupture, you must replace this part
7. Rupture disk is defective	Replace

Stopping Screen Forced through Screen Support Ring

Probable Cause	Action
1. The screen has been “unseated” from support	Carefully insert the microcarrier launch assembly into bombardment chamber to prevent movement of stopping screen. This screen movement can also occur if you close the bombardment chamber door too forcefully
2. The stopping screen support ring is damaged	Examine the support ring for evidence of scratches or dents that might prevent the screen from seating properly

Excessive Gas Usage

Probable Cause	Action
1. Leaking helium fitting in the system	Isolate the components from the He cylinder to the solenoid assembly by completely closing the metering valve (turning the knob counterclockwise until it stops). Ensure that the regulator gauge maintains pressure over time. "Soap test" the fittings by spraying a film of soapy water at the site of the fittings and looking for bubble formation. Tighten any fittings that appear to leak. Remember to close the main cylinder valve when the helium tank is not in use
2. Damaged sealing edge on gas acceleration tube (scratch or dent in metal)	Replace gas acceleration tube

Chamber Will Not Hold Vacuum

Probable Cause	Action
1. Door gasket does not seal	Test the door gasket (the large rubber O-ring around the interior perimeter of the bombardment chamber door) by spraying the outer edge of the closed door with soapy water while attempting to hold a vacuum. If the soap bubbles, replace O-ring
2. Vacuum connection at the back of the bombardment chamber is not sealed	Check for leakage (poorly fitting components) from the clamp assembly located at the rear of the unit

Sample Damage from Gas Pressure Wave

Probable Cause	Action
1. Macrocarrier not inserted	Examine and replace
2. Microcarrier launch assembly not inserted into bombardment chamber	Examine and replace
3. Incorrect sample shelf position	Reposition
4. Stopping screen absent, improperly positioned, or fails to function within the microcarrier launch assembly	Examine and replace
5. Incorrect assembly of microcarrier launch assembly components	Examine and reassemble

Unit Will Not Pressurize Gas Acceleration Tube

Probable Cause	Action
1. Rupture disk not in position or not sealed by retaining cap	Insert disk properly
2. Main cylinder valve of helium tank not open	Adjust according to Connecting the PDS-1000/He System to a Helium Source in Section 3
3. Insufficient vacuum	Examine vacuum connections (see Connecting the PDS-1000/He System to a Vacuum Source in Section 3)
4. Helium system leak	Isolate leaks in the helium pressure regulator by closing the helium cylinder valve and turning the pressure-adjusting screw one turn counterclockwise. Then note the following: <ul style="list-style-type: none"> ▪ If the high pressure gauge reading drops, there is a leak in the cylinder valve, inlet fitting, or high pressure gauge ▪ If the low pressure gauge drops, there is a leak in the valve, hose, hose fitting, outlet fitting, or low pressure gauge ▪ If the high pressure gauge drops, and at the same time the low pressure gauge rises, there is a leak in the regulator seat Defective regulators should be returned to Bio-Rad Laboratories for repair. Use of a damaged regulator, or one that has been repaired improperly is a safety hazard
5. Helium tank regulator not adjusted correctly	Adjust according to Connecting the PDS-1000/He System to a Helium Source in Section 3
6. Solenoid 3-way metering valve assembly not properly connected to main unit	Connect according to Connecting the PDS-1000/He System to a Helium Source in Section 3 and Solenoid Valve Adjustment in Section 5
7. Solenoid 3-way metering valve closed	Adjust according to Connecting the PDS-1000/He System to a Helium Source in Section 3 and Solenoid Valve Adjustment in Section 5

Section 7

Product Information

Biolistic Particle Delivery System

Catalog #	Product Description
1652256	Particle Delivery Research License
1652257	PDS-1000/He
1652258LEASE	PDS-1000 Helium Retrofit Kit , for converting the PDS-1000 to the PDS-1000/He
1652259	Voltage Converter , for countries with 220 V or 240 V
1652262	0.6 µm Gold Microcarriers , 0.25 g
1652263	1.0 µm Gold Microcarriers , 0.25 g
1652264	1.6 µm Gold Microcarriers , 0.25 g
1652278	500 Optimization Kit , includes 0.25 g each of 0.6, 1.0, and 1.6 µm gold microcarriers, 100 each of 7 rupture disks, 500 macrocarriers, 500 stopping screens
1652322	Macrocarrier Holders , set of 5
1652326	450 psi Rupture Disks , 100
1652327	650 psi Rupture Disks , 100
1652328	900 psi Rupture Disks , 100
1652329	1,100 psi Rupture Disks , 100
1652330	1,350 psi Rupture Disks , 100
1652331	1,550 psi Rupture Disks , 100
1652332	1,800 psi Rupture Disks , 100
1652335	Macrocarriers , 500
1652336	Stopping Screens , 500

Spare Parts for Helium Biolistic System Assembly (PDS-1000/He)

Spare Part #	Product Description
Plastic Case Version	
1009110	Macrocarrier Cover Lid
1652322	Microcarriers , 5
800-007	Torque Lever Assembly
8003013	Solenoid Valve Cord Assembly
8003018	Gas Acceleration Tube Assembly
8003030	Solenoid Valve
8003036	2.5' PEEK Tubing Assembly
8003039	6' PEEK Tubing Assembly
9005156	Vacuum Switch Replacement Kit
9100074	Acceleration Tube O-Ring
9100757	Retaining Spring
9101201	Set Screw
9104192	Flow Control Needle Valve
9109500	Cylinder Pressure Regulator
9109511	Helium Pressure Gauge
9109515	Vacuum Gauge
9109529	Wrench, open end, 1/4 x 5/16"
Metal Case Version (Discontinued)	
8003014	Complete Door Assembly
9109549	Complete Filter Housing
9109526	Filter Housing O-Ring

Spare Part #	Product Description
Gunpowder Version (Discontinued)	
9109538	Accelerators, 100
9203009	Locking Plug
9100019	Perimeter Seal Door O-Ring
9203061	Pin Loader
9109522	Tornado Brush
9109530	Cleaning Rod
9109531	Dud Remover Rod
9203022	Large Rubber Stopping Washer
9203023	Steel Blue Stopping Disk
9203026	Small Rubber Stopping Washer
9203035	Latch, Door 2
9203058	Mid Seal Door O-Ring
9203007	Rupture Disk Retainer Cap
9203008	Microcarrier Launch Shelf
9203010	Adjustable Nest
9203011	Fixed Nest
9203013	Stopping Screen Holder
9203014	5 mm Spacer Ring
9203015	Microcarrier Cover Lid
9203016	1/8" Gap Adjustment Tool
9203017	1/4" Gap Adjustment Tool
9203018	3/8" Gap Adjustment Tool
9203060	Petri Dish Holder
9203068	18" Vacuum Hose
9203085	Red Seals, 500

Many spare parts, accessories, and consumables for this system are no longer being manufactured. Please call your local Bio-Rad office with inquiries regarding products related to your PDS-1000/He System. In the U.S. call 1-800-4BIORAD (1-800-424-6723).

Section 8 Appendices

Cleaning the PDS-1000/He Device

Do not autoclave or flame sterilize the PDS-1000/He Instrument. Autoclaving will wet electrical components, thus creating a potential electrical shock hazard.

1. Chamber

Clean the chamber with 70% ethanol. Allow time for drying.

2. Gas Acceleration tube

The acceleration tube can be sterilized by partially immersing in 70% ethanol followed by drying in a sterile environment. Do not autoclave the gas acceleration tube. Do not immerse the pressure gauge in the sterilizing solution.

- a. Unscrew the rupture disk retaining cap from the gas acceleration tube.
- b. Disconnect the tubing assembly from the gas inlet on the rear of the acceleration tube (See Figure 1.2). Remove the screws from the hold-down brackets on the top of the instrument; two sets are located on either side of the acceleration tube. CAREFULLY remove the tube from the top of the sample chamber.

Metal Case Version

The metal chamber design has been offered by Bio-Rad Laboratories since the product line was acquired from E.I du Pont de Nemours in 1991. The original metal case design is capable of operating either the gunpowder acceleration mechanism or the helium acceleration mechanism (Figure 8.1). The gunpowder acceleration delivery system has been discontinued. The current plastic case version of the PDS-1000/He System was introduced in March 1995 and operates only with helium. The target distances and internal dimensions of the plastic case are equivalent to those found in the metal case when used with helium acceleration mechanism.



Fig. 8.1. Front view of metal case version of PDS-1000/He Instrument.

Aspects Unique to the Metal Chamber Design

- Metal bombardment chamber
- Top slots of the metal case. These are used only for the target shelf placement when using the gunpowder acceleration system
- Door construction:
Gasket design changed. Portions dedicated to use with gunpowder acceleration were removed.
Door brace added in plastic case design for additional safety
- External vacuum filter assembly components. Located on rear and side of metal case unit. These components are used to capture spent gunpowder in thin oil film inside trap after bombardment. See rear view (Figure 8.2) for identification of components
- Helium tubing is stainless steel in the metal case version and is composed of PEEK Plastic in the plastic case version
- Power cord is permanently attached in the metal case design but is detachable in the plastic case design

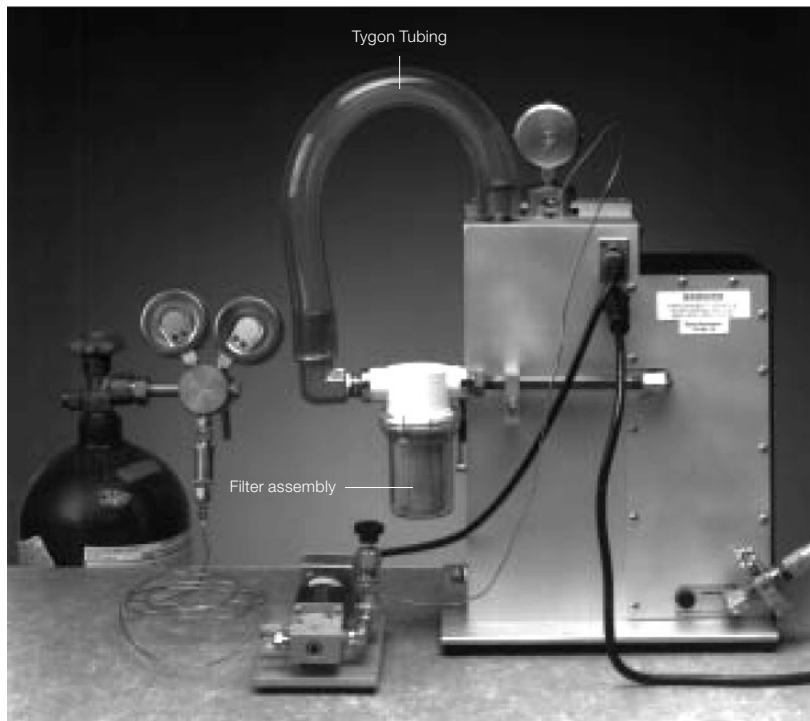


Fig. 8.2. Rear view of metal case version of PDS-1000/He System.

Attachment/Removal of Vacuum Filter Assembly for the Metal Case PDS-1000/He System

Wrap PTFE tape on the threaded brass vacuum filter fitting on the rear of the PDS-1000/He System and attach the filter assembly (Figure 8.2). The vacuum filter assembly includes a plastic filter housing and brass filter hose fitting, plus the Tygon Tubing. Turn the filter assembly approximately four complete turns, ending with the filter housing parallel to the unit and the Tygon Tubing facing upward.

Push the open end of the Tygon Tubing onto the baffle located on top of the PDS-1000/He System.

Note: The filter assembly must be firmly supported as the Tygon Tubing is being pushed onto the baffle on top of the unit (or onto the brass filter hose fitting) to avoid cracking the plastic filter housing.

Table 8.1 Sterilization and decontamination compatibility.

Method	Material						
	Aluminum	Buna-N	Delrin	Neoprene	Nylon	Polycarbonate	Polyethylene
Sterilization							
Autoclave, 121°C	U	M	M	U	U	M	M
UV irradiation	S	*	S	*	M	U	U
Ethylene oxide	S	U	S	*	S	S	S
Formaldehyde, gas	S	S	S	S	S	S	S
Gluteraldehyde, 2%	S	S	S	S	S	S	S
Biological Disinfection							
Alcohol, 70%	S	S	M	S	S	M	S
Hypochlorite, 5%	U	M	U	M	S	S	S
Gluteraldehyde, 2%	S	S	S	S	S	S	S
Phenolic derivatives	U	U	U	U	U	U	S
Formaldehyde	M	M	S	S	S	S	S

Key: S = Satisfactory; M = Moderate attack, may be satisfactory, suggest testing; U = Unsatisfactory, not recommended;
 * Performance unknown, suggest testing.

Method	Material					
	Polyvinyl Chloride	PTFE	Stainless Steel	Tygon	Acrylic	Painted Surfaces
Sterilization						
Autoclave, 121°C	M	S	S	M	U	U
UV irradiation	S	S	S	S	U	S
Ethylene oxide	S	S	S	S	S	S
Formaldehyde, gas	S	S	S	S	S	S
Gluteraldehyde, 2%	S	S	S	S	S	S
Biological Disinfection						
Alcohol, 70%	M	S	M	M	M	S
Hypochlorite, 5%	U	S	M	M	S	S
Gluteraldehyde, 2%	S	S	S	S	S	S
Phenolic derivatives	U	S	U	U	U	S
Formaldehyde	M	S	S	M	S	S

Key: S = Satisfactory; M = Moderate attack, may be satisfactory, suggest testing; U = Unsatisfactory, not recommended;
 * Performance unknown, suggest testing.

Specifications

General Specifications, PDS-1000/He Biolistic Particle Delivery System

Dimensions (W x D x H)	29 x 25.5 x 47.5 cm
Construction	Aluminum, ABS plastic, and acrylic chassis
Weight	15 kg
Electrical input voltage	100–120 VAC, 50–60 Hz
Maximum current	<5 A
Mechanical	
Fuse	6.3 A, 250 V, 5 x 20 mm
Vacuum	<0.4" Hg/min leakage
Over-pressure	0.5 psi relief valve, self-resetting
Environmental	
Operating	0–35°C (32–95°F) temperature, 0–95% noncondensing humidity
Storage	0–70°C (32–158°F) temperature, 0–95% noncondensing humidity

Source Specifications, PDS-1000/He Biolistic Particle Delivery System

The following are the sole source specifications for the Biolistic PDS-1000/He System:

1. A system capable of generating high-pressure gas delivery with a mechanism for delivering an instantaneous cold gas shock wave into an enclosure where the gas shock is released, contained, and vented.
2. The unit contains a throat region in the high pressure gas delivery system that allows for interchangeable inserts (rupture disks), which translates the cold gas shock into acceleration of microprojectiles coated with biological molecules for delivery into diverse target cells/tissue without killing the cells and/or tissue. Seven interchangeable disks are available in 450, 650, 900, 1,100, 1,350, 1,550, and 1,800 pounds per square inch (psi) pressure.
3. The system employs a method for launching the biologically coated microcarriers from the target side of the planar surface of a flexible plastic sheet that is positioned to be accelerated by the gas shock wave such that it will move freely in the direction of the target cells until restrained by a barrier. The barrier permits the microcarriers to move toward the target cells and also effectively baffles the gas shock wave to deflect some of the force of the cold gas shock wave away from the target cells with sufficient acceleration of microprojectiles coated with biological molecules for delivery into diverse target cells/tissue without killing the cells and/or tissue.

Performing a Bombardment – Quick Guide Tear-Out

Before the Bombardment

1. Select/adjust bombardment parameters for gap distance between rupture disk retaining cap and microcarrier launch assembly. Place stopping screen support in proper position inside fixed nest of microcarrier launch assembly.
2. Check helium supply (200 psi over the rated rupture pressure).
3. Clean/sterilize
 - **Equipment** – rupture disk retaining cap, microcarrier launch assembly
 - **Consumables** – macrocarriers/macrocarrier holders
4. Wash microcarriers and resuspend in 50% glycerol.
5. Coat microcarriers with DNA and load onto sterile macrocarriers/macrocarrier holders the day of experiment.

Firing the Device

1. Plug in power cord from main unit to electrical outlet.
2. Power ON.
3. Sterilize chamber walls with 70% ethanol.
4. Load sterile rupture disk into sterile retaining cap.
5. Secure retaining cap to end of gas acceleration tube (inside, top of bombardment chamber) and tighten with torque wrench.
6. Load macrocarrier and stopping screen into microcarrier launch assembly.
7. Place microcarrier launch assembly and target cells in chamber and close door.
8. Evacuate chamber, hold vacuum at desired level (minimum 5" Hg).
9. Bombard sample: hold Fire button continuously depressed until rupture disk bursts and helium pressure gauge drops to zero.
10. Release Fire button.

After the Bombardment

1. Release vacuum from chamber.
2. Remove target cells from chamber.
3. Unload macrocarrier and stopping screen from microcarrier launch assembly.
4. Unload spent rupture disk.
5. Remove helium pressure from the system (after all experiments for the day have been completed).

Go to [bio-rad.com/PDS1000/training](https://www.bio-rad.com/PDS1000/training) for additional PDS-1000/He System training and overview information, including instrument setup, consumable preparation, and instrument operation and shutdown.

Section 9

References

- Gordon-Kamm, et al. (1990). Transformation of maize cells and regeneration of fertile transgenic plants. *Plant Cell* 1990 2, 603–618.
- Sanford et al. (1993). Optimizing the biolistic process for different biological applications. *Methods Enzymol* 217, 482–509.

Visit bio-rad.com/PDS1000 for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00
Brazil 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00
Finland 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190
India 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670
The Netherlands 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717
Russia 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00
Switzerland 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

