

Bio-Plex Pro Human Diabetes Assays

Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Human Diabetes Assays	10000094511

This guide can be used to prepare and run a full 1 x 96-well assay plate. For more information on a given step, refer to the complete instruction manual. New users can go to bio-rad.com/bio-plex and download the manual, which includes detailed instructions and a list of kit components.

IMPORTANT! Pay close attention to **vortexing**, **shaking**, and **incubation** instructions. Deviation from the protocol may result in low assay signal and assay variability.

Initial Preparation

1. Plan the plate layout.
2. Start up/warm up the Bio-Plex Multiplex Immunoassay System (**30 min**).
 - Bring diluents, including wash buffer, assay buffer, standard diluent HB, detection antibody diluent HB, and sample diluent HB, to room temperature (RT). Keep the other items on ice until needed
 - Mix by inversion to ensure all salts are in solution
 - Prepare 1x wash buffer: dilute **1 part** 10x wash buffer (60 ml) with **9 parts** distilled water (540 ml)
 - Begin to thaw frozen samples
3. After thawing samples, prepare as shown.

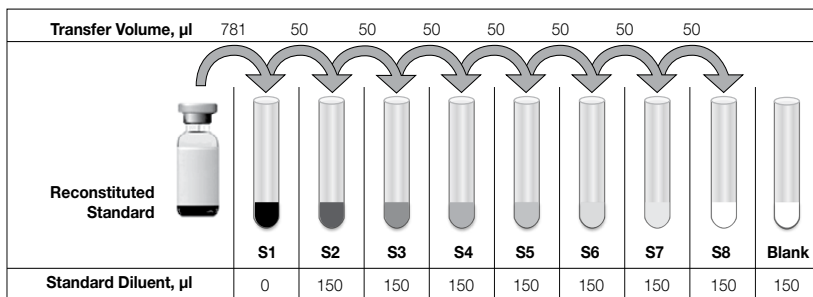
Sample Type	Diluent	Add Bovine Serum Albumin (BSA)	Sample Dilution
Serum and plasma	Sample diluent	None	Fourfold (1:4)
Culture media, with serum	Culture media	None	Neat to 1:10
Culture media, serum-free	Culture media	To 0.5% final (w/v)	Neat to 1:10

Bio-Plex Pro Human Diabetes Assays Quick Guide

Sample Type	Diluent	Add BSA	Sample Dilution
For Adiponectin Assay			
Serum and plasma	Serum-based diluent	None	Human (1:2,500)
For Adipsin Assay			
Serum and plasma	Serum-based diluent	None	Human (1:2,500)

Note: User-defined validation is required for the use of other dilution factors.

- Calibrate the Bio-Plex System by following the prompts in Bio-Plex Manager Software. This can be done now or during an assay incubation step.
- Reconstitute a single vial of standards in **781 μ l** of standard diluent HB or a diluent similar to the final sample type or matrix. **Vortex** for **5 sec** and incubate **on ice** for **30 min**.
- Prepare a fourfold standard dilution series and blank as shown. **Vortex** for **5 sec** between liquid transfers.



- Vortex** the coupled beads at medium speed for **30 sec** and **dilute to 1x** in Bio-Plex Assay Buffer as shown. Protect from light.

Premixed Panels

Number of Wells	10x Beads, μ l	Assay Buffer, μ l	Total Volume, μ l
96	570	5,130	5,700

Singleplex Assays

Number of Wells	Singleplex #1	Singleplex #2	Assay Buffer, μ l	Total Volume, μ l
	20x Beads, μ l	20x Beads, μ l		
96	285	285	5,130	5,700

Note: 20x singleplex beads allow multiplexing up to 20 analytes.

Running the Assay

Note: Make sure all assay components are at RT before proceeding.

1. **Vortex** the diluted (1x) beads for a minimum of **10 sec**. Add **50 µl** to each well of the assay plate.
2. **Wash the plate two times** with **100 µl** Bio-Plex Wash Buffer.
3. **Vortex** samples, standards, and blank. Add **50 µl** to each well.
4. Cover the plate with sealing tape and protect from light with aluminum foil. Incubate on shaker at **850 ± 50 rpm** for **1 hr** at RT.
5. With 10 min left in the incubation, **vortex** the detection antibodies for **5 sec** and quick-spin to collect liquid. **Dilute to 1x** as shown.

Premixed Panels

Number of Wells	10x Detection Antibodies, µl	10x Detection Antibody Diluent HB, µl	Total Volume, µl
96	300	2,700	3,000

Singleplex Assays

Number of Wells	Singleplex #1	Singleplex #2	Detection Antibody Diluent HB, µl	Total Volume, µl
	20x Detection Antibodies, µl	20x Detection Antibodies, µl		
96	150	150	2,700	3,000

Note: 20x singleplex beads allow multiplexing up to 20 analytes.

6. **Wash the plate three times** with **100 µl** wash buffer.
7. **Vortex** the diluted (1x) detection antibodies. Add **25 µl** to each well.
8. Cover and incubate at **850 ± 50 rpm**, as in step 4, for **30 min** at RT. Meanwhile, prepare the Bio-Plex Manager Software protocol; enter standard S1 values provided in the assay kit.
9. With 10 min left in the incubation, **vortex** the 100x streptavidin-phycoerythrin (SA-PE) for **5 sec** and quick-spin to collect liquid. **Dilute to 1x** as shown and protect from light.

Number of Wells	100x SA-PE, µl	Assay Buffer, µl	Total Volume, µl
96	60	5,940	6,000

10. **Wash the plate three times** with **100 µl** wash buffer.
11. **Vortex** the diluted (1x) SA-PE. Add **50 µl** to each well.

Bio-Plex Pro Human Diabetes Assays Quick Guide

- Cover and incubate at 850 ± 50 rpm, as in step 4, for **10 min** at RT.
- Wash the plate three times** with **100 μ l** wash buffer.
- Resuspend the beads in **125 μ l** assay buffer. Cover the plate as in step 4 and shake at 850 ± 50 rpm for **30 sec**.
- Remove the sealing tape and **read plate** using the following settings:

Instrument	RP1 (PMT)	DD Gates	Bead Events
Bio-Plex 200*	Low	5,000 (low), 25,000 (high)	50
Bio-Plex 3D*	Enhanced	Select MagPlex Beads	50
Bio-Plex MAGPIX*	N/A, use default instrument settings	N/A, use default instrument settings	Default

* Or similar Luminex System.

Bio-Rad and Bio-Plex are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

Luminex is a trademark of Luminex Corporation.

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

BIO-RAD

**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

