

Bio-Rad transforms electroporation

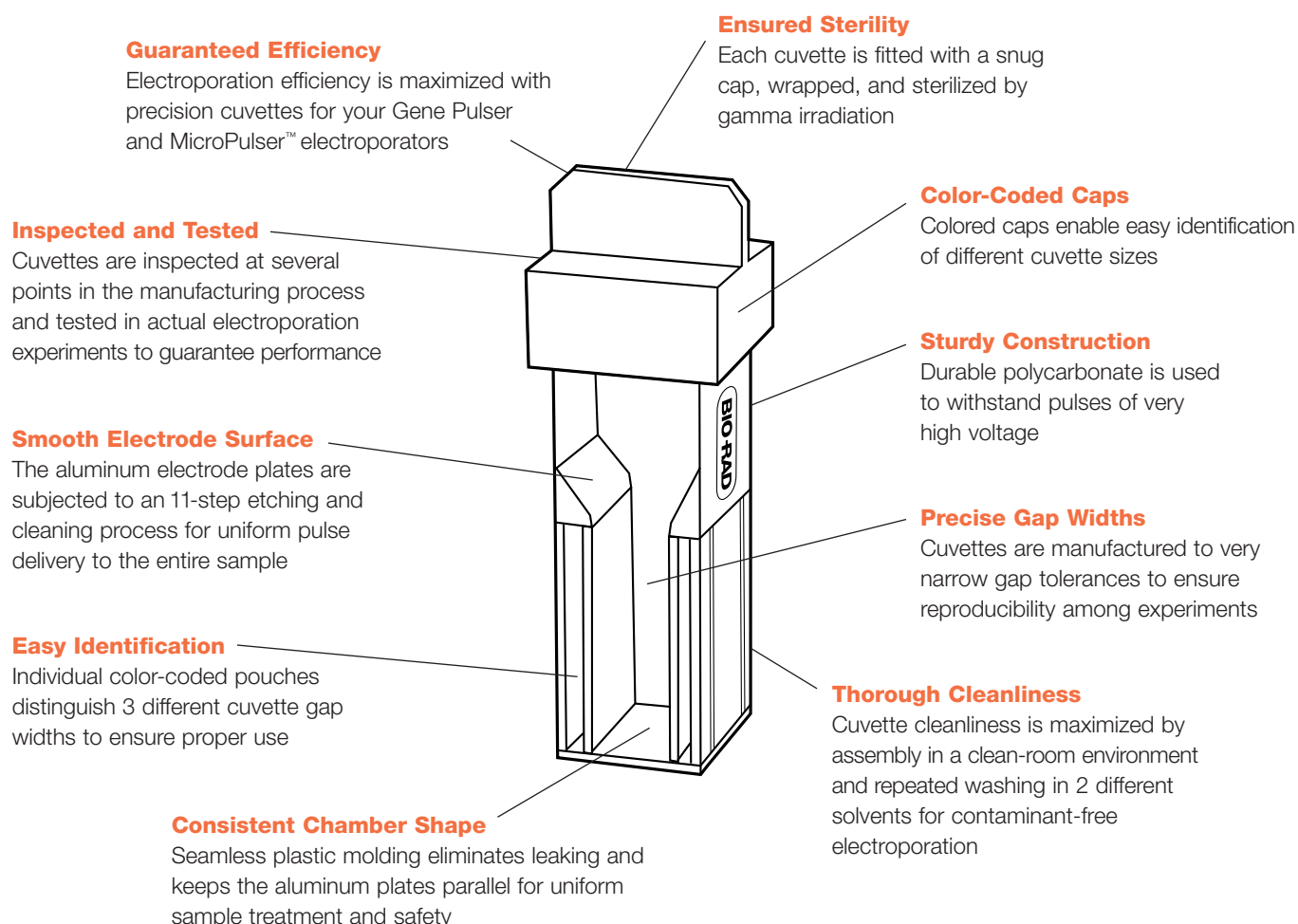


Gene Transfer Is Evolving with Gene Pulser® Cuvettes
to Maximize Electroporation Efficiencies

BIO-RAD

Bio-Rad Electroporation Cuvettes

features and benefits



Ordering Information

Catalog #	Description
165-2083	MicroPulser/Gene Pulser Cuvettes, 0.1 cm gap, 5 (mini pack)
165-2089	MicroPulser/Gene Pulser Cuvettes, 0.1 cm gap, 50 (standard pack)
165-2093	MicroPulser/Gene Pulser Cuvettes, 0.1 cm gap, 500 (jumbo pack)
165-2082	MicroPulser/Gene Pulser Cuvettes, 0.2 cm gap, 5 (mini pack)
165-2086	MicroPulser/Gene Pulser Cuvettes, 0.2 cm gap, 50 (standard pack)
165-2092	MicroPulser/Gene Pulser Cuvettes, 0.2 cm gap, 500 (jumbo pack)
165-2081	MicroPulser/Gene Pulser Cuvettes, 0.4 cm gap, 5 (mini pack)
165-2088	MicroPulser/Gene Pulser Cuvettes, 0.4 cm gap, 50 (standard pack)
165-2091	MicroPulser/Gene Pulser Cuvettes, 0.4 cm gap, 500 (jumbo pack)



**Bio-Rad
Laboratories**

Life Science
Group

Web site www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 507 6191 **Canada** (905) 712-2771 **China** 86-10-8201-1366/68 **Denmark** 45 44 52-1000 **Finland** 358 (0)9 804 2200 **France** 01 47 95 69 65 **Germany** 089 318 84-177 **Hong Kong** 852-2789-3300 **India** (91-124) 6398112/113/114 **Israel** 03 951 4124 **Italy** 34 91 590 5200 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 52 5 534 2552 to 54 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280 **Norway** 47-23-38-41-30 **Russia** 7 095 979 98 00 **Singapore** 65-2729877 **Spain** 34-91-590-5200 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 061-717-9555 **United Kingdom** 0800-181134