

# Bio-Rad transforms electroporation

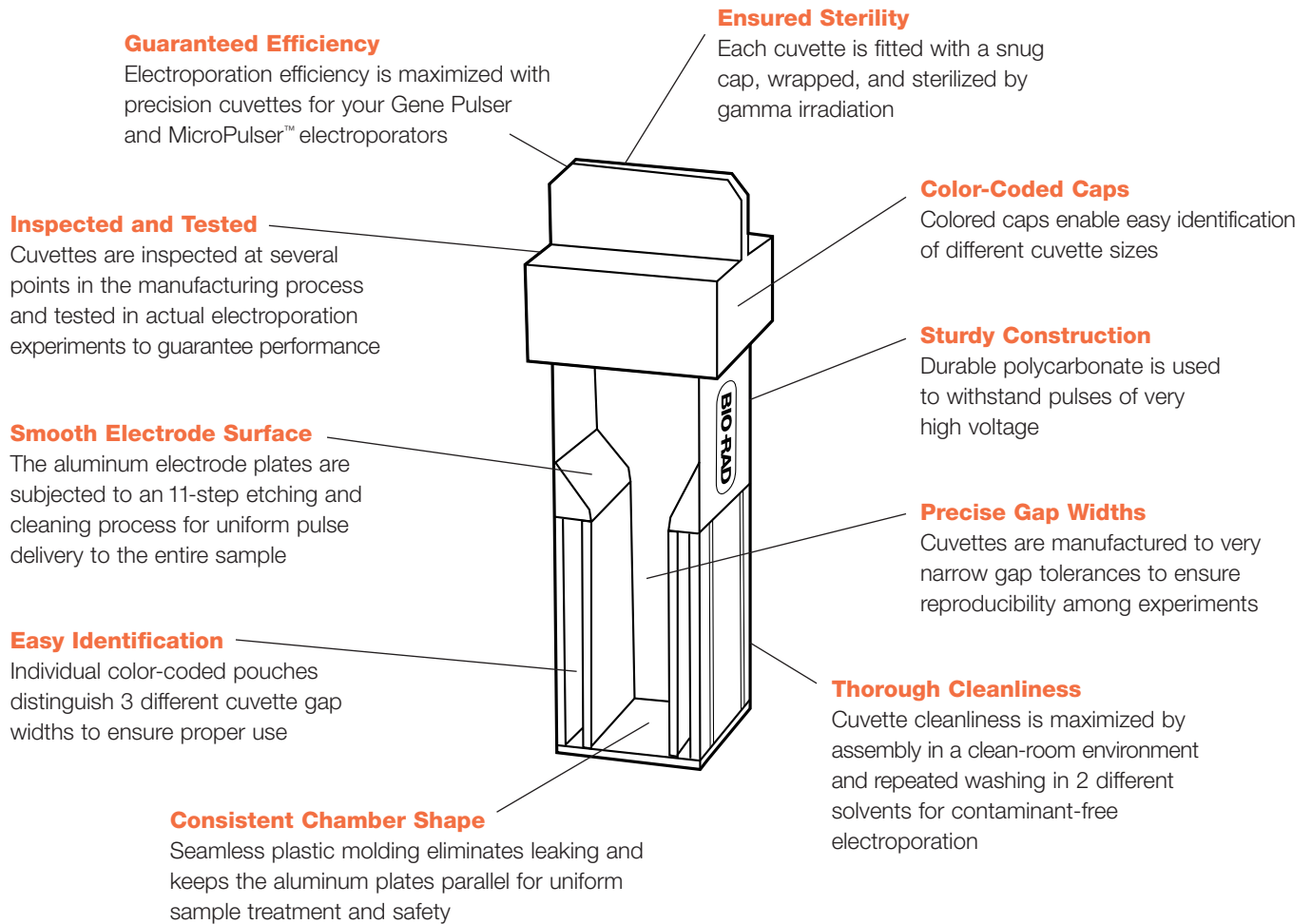


Gene Transfer Is Evolving with Gene Pulser® Cuvettes  
to Maximize Electroporation Efficiencies

**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**

**Bio-Rad  
Laboratories**

Life Science  
Group

Web site [www.bio-rad.com](http://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