

GENE TRANSFER: ELECTROPORATION

MicroPulser™ Electroporator

- Compact, space-saving design
- Audible and visible pulse indicators
- Display of time constant and actual voltage delivered to monitor reproducibility
- 3,000 V capability for improved efficiency in cuvettes with larger volume

An Ideal Gene Transfer Tool for Bacteria, Fungi, and Other Microorganisms

Advantages of Electroporation

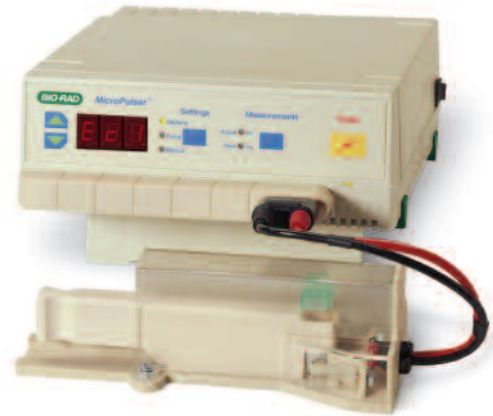
The MicroPulser electroporator is a simple yet versatile electroporator that provides a safe and reproducible way for you to transform bacteria, fungi, and other microorganisms. Transformation efficiencies much higher than those obtained by chemical methods can be achieved by electroporation.

The MicroPulser electroporator's compact, space-saving design comes with preset optimized programs for transformation of commonly studied bacteria and fungi. Its audible and visible pulse indicators display time constant and actual voltage delivered (200–3,000 V with 10 V precision) to monitor reproducibility and ensure accuracy, while its arc-quenching (ARQ) system protects valuable samples. A broad range of parameters is available for manual optimization.

The MicroPulser electroporator offers safe operation, rapid optimization, and outstanding results.

Unique Features

- ARQ system that reduces arcing, protecting against loss of valuable samples
- Rapid program selection — preset optimized programs for commonly studied bacteria and fungi
- Efficient sample handling with one-button pulse delivery, rapid charge time, and attached cuvette chamber



BIO-RAD

MicroPulser Electroporator

Preprogrammed Settings

The MicroPulser is preprogrammed with settings for electroporation of a number of commonly used organisms. The table below lists programs that are available under Bacteria Settings and Fungi Settings.

Mnemonic	Organism	Parameters		
		Voltage (kV)	Number of Pulses	Time Constant* (ms)
Bacteria				
Ec1	<i>E. coli</i> (0.1 cm cuvette)	1.8	1	—
Ec2	<i>E. coli</i> (0.2 cm cuvette)	2.5	1	—
StA	<i>S. aureus</i> (0.2 cm cuvette)	1.8	1	2.5
Agr	<i>A. tumefaciens</i> (0.1 cm cuvette)	2.2	1	—
Ec3	<i>E. coli</i> (0.2 cm cuvette)	3.0	1	—
Fungi				
Sc2	<i>S. cerevisiae</i> (0.2 cm cuvette)	1.5	1	—
Sc4	<i>S. cerevisiae</i> (0.2 cm cuvette)	3.0	1	—
ShS	<i>S. pombe</i> (0.2 cm cuvette)	2.0	1	—
Dic	<i>D. discoideum</i> (0.4 cm cuvette)	1.0	2	1.0
Pic	<i>P. pastoris</i> (0.2 cm cuvette)	2.0	1	—

* Unless the pulse time is truncated below 5 ms, the unit will deliver the optimal time constant of ~5 ms to samples in high-resistance media.

Cuvettes

Bio-Rad's high-quality electroporation cuvettes are available in three different gap widths — 0.4, 0.2, and 0.1 cm — for optimal field strength delivery to a wide range of cell types. The cuvettes ensure maximum electroporation efficiency and reproducibility between experiments, are color-coded, and have been sterilized by gamma irradiation.

Specifications

Input voltage	100–120 V or 220–240 V
Maximum output voltage and current	3,000 V peak into >600 W load; limited at 100 A peak maximum
Output waveform	Decaying or truncated-decaying exponential waveform with RC time constant
Output voltage and pulse duration adjustment	Voltage adjustable in 200–3,000 V range with 10 V precision; 5 ms default or 1–4 ms with 0.1 ms precision; 5 bacterial and 5 fungal preset programs
Ambient operating temperature	3.5–35°C
Dimensions (W x D x H)	29 x 21 x 8 cm
Weight	2.9 kg

Ordering Information

Catalog # Description

MicroPulser Electroporator

165-2100 **MicroPulser Electroporator**, includes a cuvette chamber with leads, 10 sterile cuvettes (5 each of 0.1 cm and 0.2 cm gap)

Cuvettes

Standard Packs

165-2088 Gene Pulser®/MicroPulser Cuvettes, 0.4 cm gap, 50 (regular pack)
 165-2086 Gene Pulser/MicroPulser Cuvettes, 0.2 cm gap, 50 (regular pack)
 165-2089 Gene Pulser/MicroPulser Cuvettes, 0.1 cm gap, 50 (regular pack)

Jumbo Packs*

165-2091 Gene Pulser/MicroPulser Cuvettes, 0.4 cm gap, 500 (jumbo pack)
 165-2092 Gene Pulser/MicroPulser Cuvettes, 0.2 cm gap, 500 (jumbo pack)
 165-2093 Gene Pulser/MicroPulser Cuvettes, 0.1 cm gap, 500 (jumbo pack)

Mini Packs

165-2081 Gene Pulser/MicroPulser Cuvettes, 0.4 cm gap, 5 (mini pack)
 165-2082 Gene Pulser/MicroPulser Cuvettes, 0.2 cm gap, 5 (mini pack)
 165-2083 Gene Pulser/MicroPulser Cuvettes, 0.1 cm gap, 5 (mini pack)

* Please inquire about volume pricing.

For more information, visit us on the Web at
www.bio-rad.com/genetransfer/ or
www.bio-rad.com/rnai/

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

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 3237 9400
 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050
 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280
 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723
 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000