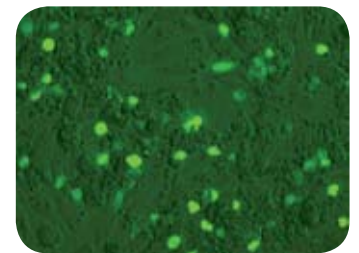


Primary Transfection
Tool

Gene Pulser MXcell™ Electroporation System

Obtain better gene delivery while maintaining cell viability in mammalian cells, especially difficult-to-transfect cells such as primary and stem cells.

- Preprogrammed protocols to quickly determine optimal starting conditions, regardless of cell type
- Manual programming of protocols to account for the wide variations in primary cells and stem cells
- Choice of delivery format — cuvette or plate — to electroporate a limited number of primary cells or a larger number when scaling up
- 1–24 different protocols on a single electroporation plate — control the amount of cell culture required to obtain sufficient cell numbers for analysis
- Fast pulse time to minimize handling of cells
- Compatibility with Gene Pulser® electroporation buffer to ensure efficient gene delivery, while maintaining cell viability



Mouse embryonic stem (ES) cells electroporated using the Gene Pulser MXcell system and Gene Pulser electroporation buffer.

Postelectroporation mouse ES cells were grown on STO feeder cells for visualization by fluorescent microscopy.

BIO-RAD

The Primary Solution for Your Transfection Needs

The MXcell is an affordable electroporation system offering a choice of delivery tools. The system consists of a power module to generate a pulse, a plate chamber, and a 96-well electroporation plate. The optional Gene Pulser MXcell™ ShockPod™ cuvette chamber allows an easy transition between electroporation plates and cuvettes.

Flexible Electroporation System

- Preset and gradient protocols reduce programming
- Cross-compatible protocols allow a simple transition between plates and cuvettes
- Exponential and square waveforms ensure optimal results
- Choose a different parameter for each well set — up to 24 different protocols on a single electroporation plate

Multiple Delivery Options

- 96-well electroporation plate for small numbers of cells
- 24- and 12-well electroporation plates for larger laboratory-scale gene expression experiments
- Cuvettes for single transfections
- Highly reproducible results, regardless of delivery format

Format	Well Volume	Cell Concentration	Number of Well Sets
96-well plate	100–200 µl	1 x 10 ⁵ to 2 x 10 ⁶	24
24-well plate	500–800 µl	5 x 10 ⁵ to 8 x 10 ⁶	24
12-well plate	1.0–1.5 ml	1 x 10 ⁶ to 1.5 x 10 ⁷	12
0.4 cm cuvette	400–800 µl	4 x 10 ⁵ to 8 x 10 ⁶	N/A

Self-Contained Unit

- No computer or power adaptor required
- User-friendly digital interface with easy, intuitive programming controls all parameters

Specifications

Outputs

Waveform	Exponential-decay or square-wave
Voltage	10–500 V
Capacitance	10–500 V, 25–2,475 µF in 25 µF increments
Resistance (parallel)	50–1,500 Ω in 50 Ω increments, plus infinity
Sample resistance	10 Ω minimum at 10–500 V
Square-wave timing	10–500 V: 0.05–9,999.95 ms pulse duration, 1–3 pulses; 0.1–10 sec pulse interval

General

Input voltage	100–120 VAC or 220–240 VAC, 50 or 60 Hz
Power	Maximum 240 W (during short charging periods)
Operating environment	Temperature 0–35°C; humidity 0–95% (noncondensing)
Regulatory	Safety EN 61010, EMC EN 61326 Class A
Dimensions (W x D x H)*	31 x 30 x 14 cm
Weight	6.62 kg

*Includes power module and chamber.

Ordering Information

Catalog #	Description
165-2670	Gene Pulser MXcell Electroporation System , 100–240 V, 50/60 Hz, exponential-decay and square-wave delivery, includes power module, plate chamber, 1 x 96-well electroporation plate, instructions
165-2674	Gene Pulser MXcell Electroporation System With MXcell ShockPod , 100–240 V, 50/60 Hz, exponential-decay and square-wave delivery, includes power module, plate chamber, 1 x 96-well electroporation plate, Gene Pulser MXcell ShockPod cuvette chamber, instructions
165-2671	Power Module
165-2672	Plate Chamber
165-2681	96-Well Electroporation Plate
165-2682	24-Well Electroporation Plate
165-2683	12-Well Electroporation Plate
165-2673	Gene Pulser MXcell ShockPod Cuvette Chamber
165-2081	Gene Pulser/MicroPulser™ Cuvettes , 0.4 cm gap, 5
165-2088	Gene Pulser/MicroPulser Cuvettes , 0.4 cm gap, 50
165-2091	Gene Pulser/MicroPulser Cuvettes , 0.4 cm gap, 500
165-2676	Gene Pulser Electroporation Buffer , 10 x 1.8 ml
165-2677	Gene Pulser Electroporation Buffer , 30 ml

For more product information, please visit us on the Web at www.bio-rad.com/genetransfer/mxcell/.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 364 3435 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 Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000