



We Make It Easy

EasyShock Electro-Competent Cells



Simplify your research with EasyShock10B electro-competent cells, provided in easy-to-use prefilled cuvettes.

BIO-RAD



Convenience that simplifies your research

Introducing the EasyShock Format

Bio-Rad combines the high performance and versatility of its EP-Max™10B electro-competent cells with high-quality, precision-built 0.1 cm gap electroporation cuvettes to provide the ultimate convenience in electrotransformation. Achieve transformation efficiencies of $>1 \times 10^{10}$ CFU/ μ g DNA with no microtubes and only one pipetting step.

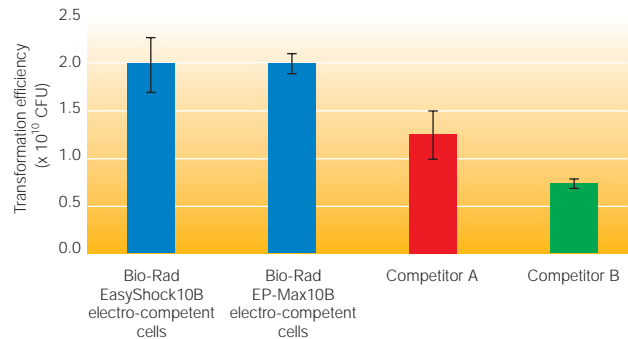
- **Easy protocol** — Simply add DNA and electroporate
- **Higher throughput** — Perform twice the transformations in half the time of a standard protocol
- **Excellent efficiency** — Achieve transformation efficiencies of $>1 \times 10^{10}$ CFU/ μ g DNA
- **Lower cost** — Even with cuvettes included, the EasyShock format is a better value than other “convenience” formats

Comparison of EasyShock and Standard Protocols

EasyShock Protocol	Standard Protocol
1. Thaw EasyShock cuvette on ice.	1. Chill cuvette and microtube on ice.
2. Add DNA to cells in cuvette and mix.	2. Thaw cells on ice.
3. Electroporate.	3. Transfer cells to chilled microtube.
4. Add SOC medium and plate cells.	4. Add DNA to cells and mix.
	5. Transfer suspension to cuvette.
	6. Electroporate.
	7. Add SOC medium and plate cells.

EasyShock Components

Cell line	EP-Max10B
Genotype	<i>F⁻ mcrA Δ(mrr-hsdRMS-mcrBC) φ80dlacZΔM15 ΔlacX74 deoR recA1 endA1 araD139 Δ(ara, leu)7697 galU galK rpsL nupG λ⁻</i>
Volume	20 μ l per cuvette
Cuvette	0.1 cm gap



A comparison of transformation efficiencies between Bio-Rad EasyShock10B, Bio-Rad EP-Max10B, and two competitors' products. Electrotransformations were performed using pUC19 control plasmid and a MicroPulser™ electroporator according to the manufacturer's recommendations.

Ordering Information

Catalog #	Description
Electro-Competent Cells — EasyShock Format	
170-3333	EasyShock10B Electro-Competent Cells, 0.1 cm electroporation cuvettes prefilled with 20 μ l of EP-Max10B cells, enough for 10 transformations
Electro-Competent Cells — Standard Format	
170-3330	EP-Max10B Electro-Competent Cells, 5 x 0.10 ml
Empty Electroporation Cuvettes	
165-2083	Gene Pulser®/MicroPulser Cuvettes, 0.1 cm gap, 5 (mini pack)
165-2089	Gene Pulser/MicroPulser Cuvettes, 0.1 cm gap, 50 (regular pack)
165-2093	Gene Pulser/MicroPulser Cuvettes, 0.1 cm gap, 500 (jumbo pack)
Electroporation Instruments	
165-2100	MicroPulser Electroporator
165-2662	Gene Pulser Xcell™ Microbial System



**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 2527 3454 **Canada** (905) 712-2771 **China** (86-21) 63052255 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84-0 **Hong Kong** 852-2789-3300 **Hungary** 36 1 455 8800 **India** (91-124)-6398112/113/114, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 55-52-00-05-20 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 331 99 99 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-6415 3188 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000