



Gene Pulser® Electroprotocols

* We recommend adapting this protocol to use the Gene Pulser electroporation buffer (catalog #165-2676, 165-2677), which increases cell viability and transfection efficiency in mammalian cell lines.

Cell type: Mouse embryonic stem cells
Strain: 129/OLA, subclone E14
Molecules: Linearized and circular DNA

Before the Pulse

Cell growth

medium: Complete GMEM consists of 1x GMEM medium (Sigma, #G5154) supplemented with 2 mM glutamine, 1 mM sodium pyruvate, 50 μ M β -mercaptoethanol, 1x nonessential amino acids, 10% (v/v) fetal bovine serum, and 500–1,000 U/ml of leukocyte inhibitory factor (LIF; Chemicon).

Wash solution: PBS
Growth phase at harvest: 90% confluent
Pre-pulse incubation: None

The Pulse

Electroporation temperature: Ambient
Electroporation medium*: PBS
Cell density: 4×10^6 to 6×10^7 cells
Volume of cells: 0.7 ml
DNA: 5.0–125 μ g
DNA resuspension buffer: PBS
Volume of DNA: 100 μ l
Instruments used: Gene Pulser Xcell with CE module

Cuvette gap: 0.4 cm
Voltage: 795 V
Field strength: —
Capacitance: 10 μ F
Resistance: Infinity
Time constant: 0.2 ms

Relevant comments: electroporation carried out in Time Constant mode using settings of 800 V, 0.2 ms time constant, and 4 mm cuvette.

After the Pulse

Cells are allowed to recover for at least 20 min at ambient temperature, diluted in complete GMEM, and plated out at a density of $2\text{--}3 \times 10^6$ cells per 10 cm dish.

Selection used: G418
Electroporation efficiency: Not determined
Percent survival: About 50%

Protocol Number: 207

Author's Name: Roland Russnak
Institution: University of California, San Francisco