



Gene Pulser® Electroprotocols

Cell Type Other Cell Types
Species Used Hydra cells, *Cnidaria*

Molecules Electroported Analoge cytoskeletal proteins (like tubulin).

Before the Pulse

Cell growth medium (Dissociation buffer) : 1.2 mM MgSO₄, 6 mM CaCl₂, 3.6 mM KCl, 6 mM pyruvate; 6 mM Na-Citrate, 12.5 mM TES buffer, 6 mM glucose, pH 6.9.
Wash solution Not given

Growth phase at harvest Not given
Pre-pulse incubation Not given

The Pulse

Electroporation Temperature Room temperature
Electroporation Medium The same as cell growth medium

Cell Density 10 (5) cells / ml
Volume of Cells 10 (4) cells / pulse

DNA Concentration Not given
DNA Resuspension Buffer Not given
Volume of DNA Not given

Instruments Used Gene Pulser® apparatus
Cuvette Gap 0.4 cm
Voltage 0.230 kV
Field Strength 0.575 kV/cm
Capacitor 25 μF
Resistor (Pulse Controller) none Ω
Time Constant 4.3 msec

After the Pulse

Outgrowth Medium The same

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature Room temperature
Length of Incubation Not given
Selection Method or Assay Used These cells are "primary cell cultures"
Electroporation Efficiency Not given
Per Cent Survival Not given

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 3/19/91
Survey Number 189