



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	Mammalian, suspension	Molecules	Fluorescent dyes (SBFI and MQAE);
Species	Rat, submandibular acini (secretory	Electroporated	(3)H-inositol.
Used	cells).		

Before the Pulse

Cell growth medium	KRH + Ca	Growth phase at harvest	Not applicable
Wash solution	Cell growth medium	Pre-pulse incubation	A variation of the growth medium, but high in potassium and low in sodium and calcium.

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Not given
Electroporation Medium*	Same as pre-pulse	Cuvette Gap	0.4 cm
Cell Density	10 mg protein / ml	Voltage	0.6 to 1.25 kV
Volume of Cells	10 mg protein / ml	Field Strength	1.5 to 3.125 kV/cm
DNA Concentration	1 mg protein / ml	Capacitor	25 μ F
DNA Resuspension Buffer	Same as Pre-Pulse buffer or KRH + Ca	Resistor	(Pulse Controller) Ω none. NOT RECOMMENDED (SEE
Volume of DNA	Not assayed	Time Constant	0.4 msec

After the Pulse

Outgrowth Medium	Not applicable	Relevant Publications and/or Comments
Outgrowth Temperature	Not given	Note: exponential values designated in parentheses.
Length of Incubation	Not given	**It is NOT RECOMMENDED to use high voltage with out the Pulse Controller.
Selection Method or Assay Used	Not given	
Electroporation Efficiency	Not given	
Per Cent Survival	Not given	

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 8/22/90
Survey Number 147
© Bio-Rad Laboratories, 1993