



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	DNA: immunoglobulin genes in pSV
Species Used	Mouse, NSO, myeloma cells	Electroporated	vector

Before the Pulse

Cell growth medium	RPMI + 10% Fetal Calf Serum (GIBCO/BRL, Sigma)	Growth phase at harvest	Not given
Wash solution	Not given	Pre-pulse incubation	10 min. on ice

The Pulse

Electroporation Temperature	0 °C	Instruments Used	Not given
Electroporation Medium*	Not given	Cuvette Gap	0.2 cm
Cell Density	12 x 10 ⁽⁶⁾ / ml	Voltage	0.250 kV
Volume of Cells	400 µl	Field Strength	1.25 kV/cm
DNA Concentration	Not given	Capacitor	960 µF
DNA Resuspension Buffer	Not given	Resistor	(Pulse Controller) Ω none
Volume of DNA	20 µl	Time Constant	0.4 to 0.8 msec

After the Pulse

Outgrowth Medium	Not given	Relevant Publications and/or Comments	Note: exponential values designated in parentheses.
Outgrowth Temperature	0 °C for 10 min. (ice)		
Length of Incubation	Not given		
Selection Method or Assay Used	mycophenolic acid		
Electroporation Efficiency	None		
Per Cent Survival	40 to 50 %		

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 5/8/92
Survey Number 140
© Bio-Rad Laboratories, 1993