



Gene Pulser® Electroprotocols

Cell Type Fungal / Yeast

Species Used *Aspergillus nidulans*

Molecules Electroported DNA: integrative plasmid

Before the Pulse

Cell growth medium Not given

Growth phase at harvest Not given

Pre-pulse incubation Not given

Wash solution Not given

The Pulse

Electroporation Temperature Not given

Electroporation Medium 1.2 M Sorbitol, 7mM NaPO₄ (pH7.2), 1 mM MgSO₄

Cell Density 4 x 10⁶ (6) protoplasts / ml

Volume of Cells Not given

DNA Concentration Not given

DNA Resuspension Buffer Not given

Volume of DNA Not given

Instruments Used Gene Pulser® apparatus

Cuvette Gap 0.2 cm

Voltage 0.400 kV, 0.700 kV

Field Strength 2.0 kV/cm, 3.5 kV/cm

Capacitor 25 μF

Resistor (Pulse Controller) Ω none

Time Constant 5.2 msec / 3.3 msec

After the Pulse

Outgrowth Medium Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Outgrowth Temperature Not given

Length of Incubation Not given

Selection Method or Assay Used Not given

Electroporation Efficiency 3.5, 4.0 transformants / μg DNA

Per Cent Survival Not given

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 3/7/91

Survey Number 170

© Bio-Rad Laboratories, 1993