



Gene Pulser® Electroprotocols

Cell Type Fungal / Yeast

Species Used *Pichia pastoris* GTS115

Molecules Electroported DNA: plasmid, pHILD2 & D4, linearized, 8 to 10 kB.

Before the Pulse

Cell growth medium Yeast Extract Potato Dextrose, YEPD, (DIFCO)

Growth phase at harvest O.D. (600) = 1.3

Pre-pulse incubation 5 minutes, 4 °C

Wash solution Cold water two times; then 1 M sorbitol, one time

The Pulse

Electroporation Temperature 25 °C but sample & cuvette at 4 °C
Electroporation Medium 1 M sorbitol

Instruments Used Gene Pulser® apparatus & Pulse Controller

Cell Density 300x concentration from harvest density [O.D. (600) = 1.3, above]

Cuvette Gap 0.2 cm

Volume of Cells 50 µl

Voltage 1.5 kV

Field Strength 7.5 kV/cm

DNA Concentration 0.5 to 2 µg / pulse

Capacitor 25 µF

DNA Resuspension Buffer 1 M sorbitol

Resistor (Pulse Controller) 400 Ω

Volume of DNA 1 to 5 µl

Time Constant approximately 8.0 msec

After the Pulse

Outgrowth Medium Minimal salts plus dextrose (MD)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. This is essentially the method described for *Saccharomyces cerevisiae* by Becker and Guarente, *Methods in Enzymol.*, **194**, 182-187(1991).

Outgrowth Temperature 30 °C

Length of Incubation 3 to 5 days

Selection Method or Assay Used Complimentation of histidine auxotrophy

Electroporation Efficiency approx. 1000 transformants / µg DNA

Per Cent Survival Not tested

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 9/2/92

Survey Number 205

© Bio-Rad Laboratories, 1993