



Gene Pulser® Electroprotocols

Cell Type Fungal / Yeast
Species Used *Saccharomyces cerevisiae* - lines derived from S288C

Molecules Electroporated DNA: pUC19 and Bluescript™ -based plasmids.

Before the Pulse

Cell growth medium YEPD (ATCC#1202/1245)

Growth phase at harvest 0.5 to 1.0 at O.D.600; ~100 mls log phase culture concentrated 200x, in

Pre-pulse incubation None required

Wash solution Water

The Pulse

Electroporation Temperature Room temperature

Instruments Used Gene Pulser® apparatus & Pulse Controller

Electroporation Medium Water

Cell Density 6×10^{10} cells / ml

Cuvette Gap 0.2 cm

Volume of Cells 40 to 50 μ l

Voltage 0.6 kV

DNA Concentration Total DNA = 0.2 to 1.0 μ g

Field Strength 3.0 kV/cm

DNA Resuspension Buffer Not given

Capacitor 25 μ F

Volume of DNA $\leq 5 \mu$ l

Resistor (Pulse Controller) 200 Ω

Time Constant Not given

After the Pulse

Outgrowth Medium Add water to final vol. 150-200 μ l. Plate on selective synthetic media (SD)

Relevant Publications and/or Comments

Note: exponential values designated in parentheses. Washed and concentrated cells can be stored by adding 80% glycerol to a final concentration of 20% and freezing at -80°C. To electroporate thawed cells: pellet cells, remove glycerol containing supernatant, wash cells in 0.5-1.0 ml water, and resuspend cells in water to original volume. Efficiency is not affected significantly by freezing but you must remove glycerol.

Outgrowth Temperature 30 °C

Length of Incubation 2 days

Selection Method or Assay Used Amino acid or nucleotide dropout

Electroporation Efficiency 2 to 10,000 transfectants / μ g DNA

Comment: removing glycerol after thawing cells may help by simply removing lysed cell contents that would alter media conductivity - this could impact efficiency (alters the time constant, t).

Per Cent Survival Not given

Name of Submitter
Institution Address

Telephone Number

Fax Number

Date Submitted 9/14/90

Survey Number 179

© Bio-Rad Laboratories, 1993