

# Introduction of Proteins into Cells by Electroporation

- ED Berglund, D. L. and Starkey, J. R., Isolation of viable tumor cells following introduction of labeled antibody to an intracellular oncogene product using electroporation, *J. Immunol. Meth.*, **125**, 79-87 (1989).
- ED Chakrabarti, R., Wylie, D. E. and Schuster, S. M., Transfer of monoclonal antibodies into mammalian cells by electroporation, *J. Biol. Chem.*, **264**, 15494-15500 (1989).
- GP+CE Cortés, F. and Ortiz, T., Induction of chromosomal aberrations in the CHO mutant EM9 and its parental line AA8 by *EcoRI* restriction endonuclease: electroporation experiments, *Mutat. Res.*, **246**, 221-226 (1991).
- GP+CE Dagher, S. F., Conrad, S. E., Werner, E. A. and Patterson, R. J., Phenotypic conversion of TK-deficient cells following electroporation of functional TK enzyme, *Exp. Cell Res.*, **198**, 36-42 (1992).
- UK Elgizoli, M., Dai, Y., Kempf, C., Koblet, H. and Michel, M. R., Semliki Forest virus capsid protein acts as a pleiotropic regulator of host cellular protein synthesis, *J. Virol.*, **63**, 2921-2928 (1989).
- GP+CE Graziadei, L., Burfeind, P. and Bar-Sagi, D., Introduction of unlabeled proteins into living cells by electroporation and isolation of viable protein-loaded cells using dextran-fluorescein isothiocyanate as a marker for protein uptake, *Anal. Biochem.*, **194**, 198-203 (1991).
- GP+CE Lambert, H., Pankov, R., Gauthier, J. and Hancock, R., Electroporation-mediated uptake of proteins into mammalian cells, *Biochem. Cell Biol.*, **68**, 729-734 (1990).
- GP+CE Langlet, C. and Schmitt-Verhulst, A.-M., Electroporation of CTL clones: a useful method to investigate signalling pathways leading to the expression of effector functions, *J. Immunol. Meth.*, **151**, 107-115 (1992).
- ED Morgan, W. F., Chung, H. W., Phillips, J. W. and Winegar, R. A., Restriction endonucleases do not induce sister-chromatid exchanges in Chinese hamster ovary cells, *Mutat. Res.*, **226**, 203-209 (1989).
- ED Negrutskii, B. S. and Deutscher, M. P., Channeling of aminoacyl-tRNA for protein synthesis *in vivo*, *Proc. Natl. Acad. Sci. USA*, **88**, 4991-4995 (1991).
- ED Tsongalis, G. J., Lambert, W. C. and Lambert, M. W., Correction of the ultraviolet light induced DNA-repair defect in xeroderma pigmentosum cells by electroporation of a normal human endonuclease, *Mutat. Res.*, **244**, 257-263 (1990).
- SW Uno, I., Fukami, K., Kato, H., Takenawa, T. and Ishikawa, T., Essential role for phosphatidylinositol 4,5-bisphosphate in yeast cell proliferation, *Nature*, **333**, 188-190 (1988).
- ED Wilson, A. K., Claypool, W. D. and de Lanerolle, P., The use of electroporation (EP) to insert antibodies to myosin light chain kinase (MK-Ab) into rat alveolar macrophages (AM), *Biophys. J.*, **51**, 214a (1987).
- ED Wilson, A. K., Claypool, W. D. and de Lanerolle, P., Incorporation of affinity-purified antibodies to myosin light chain kinase (MLCK) by electroporation (EP) inhibits macrophage (M $\phi$ ) motility, *Fed. Proc.*, **46**, 1097 (1987).
- ED Winegar, R. A., Phillips, J. W., Youngblom, J. H. and Morgan, W. F., Cell electroporation is a highly efficient method for introducing restriction endonucleases into cells, *Mutat. Res.*, **225**, 49-53 (1989).
- ED Yorifuji, T. and Mikawa, H., Co-transfer of restriction endonucleases and plasmid DNA into mammalian cells by electroporation: effects on stable transformation, *Mutat. Res.*, **243**, 121-126 (1990).
- SW Zeira, M., Tosi, P.-F., Mouneimne, Y., Lazarte, J., Sneed, L., Volsky, D. J. and Nicolau, C., Full-length CD4 electroinserted in the erythrocyte membrane as a long-lived inhibitor of infection by human immunodeficiency virus, *Proc. Natl. Acad. Sci. USA*, **88**, 4409-4413 (1991).

## Key to Electroporation Units Used

- GP** Gene Pulser<sup>®</sup> apparatus
- CE** Capacitance Extender
- ED** Exponential Decay instrument (other than a Gene Pulser apparatus)
- SW** Square Wave instrument
- UK** Unknown

**BIO-RAD****Bio-Rad  
Laboratories****Life Science  
Group**

2000 Alfred Nobel Drive  
Hercules, California 94547  
Telephone (510) 741-1000  
Fax: (510) 741-1060

**Eastern Regional Office**, 85A Marcus Dr., Melville, New York 11747 • Phone (516) 756-2575 • Fax (516) 756-2594  
**European Headquarters**, Bio-Rad Laboratories, Dreve du Sénéchal, 19, B-1180 Brussels • Phone 02 375 59 70 • Fax 02 374 61 62  
**Australia**, Bio-Rad Laboratories Pty Limited, Unit 11, 112-118 Talavera Rd P.O. Box 371, North Ryde, N.S.W. 2113 • Phone 02-805-5000 • Fax 02-805-1920  
**Austria**, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, A-1130 Wien • Phone 0222-877 89 01 • Fax 0222-876 56 29  
**Belgium**, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, B-9810 Nazareth Eke • Phone 091-85 55 11 • Fax 091-85 65 54  
**Canada**, Bio-Rad Laboratories (Canada) Ltd., 5149 Bradco Boulevard, Mississauga, Ontario L4W 2A6 • Phone (416) 624-0713 • Fax (416) 624-3019  
**China**, Bio-Rad Laboratories, Yanshan Hotel Office Tower, #1307, 138A Haidian Road, Beijing • Phone 2563146 • Fax 2564308  
**France**, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94203 Ivry Sur Seine Cedex • Phone 01-49 60 68 34 • Fax 01-46 71 24 67  
**Germany**, Bio-Rad Laboratories GmbH, Heidemannstraße 164, Postfach 45 01 33, D-8000 München 45 • Phone 089-318 84-0 • Fax 089-318 84 100  
**Italy**, Bio-Rad Laboratories S.r.l., Via Cellini, 18A, 20090 Segrate Milano • Phone 02-21609.1 • Fax 02-21609-399  
**Japan**, Nippon Bio-Rad Laboratories, K. K., Sumitomo Seimei Kachidoki Bldg 5-3-6 Kachidoki, Chuo-Ku, Tokyo 104 • Phone 03-3534-7515 • Fax 03-3534-8027  
**The Netherlands**, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 08385-40666 • Fax 08385-42216  
**New Zealand**, Bio-Rad Laboratories, Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097  
**Pacific**, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257  
**Scandinavia**, Bio-Rad Laboratories, Kanalvägen 10C, 19461 Upplands Väsby, Sweden • Phone 46 (0) 8 590-73489 • Fax 46 (0) 8 590-71781  
**Spain**, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98  
**Switzerland**, Bio-Rad Laboratories AG, Kanalstrasse, 17, 8152 Glattpfugg • Phone 01-810 16 77 • Fax 01-810 19 33  
**United Kingdom**, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Phone 0800 181134 • Fax 0442 259118