

PEG: Polyethylene Glycol Solution

CATALOG#: ENZ-70007-0001

1 ml/vial

PRODUCT DESCRIPTION

This product is a 50% (w/v), average mol wt 1,450, prescreened and sterile-filtered solution in DPBS without calcium. PEG induces hybridization and promotes fusion of the cell membranes of B cells with myeloma fusion partner.

STABILITY AND STORAGE

Product stable at 2-8°C until expiration date as indicated on label. Product is sterility tested.

DIRECTIONS FOR USE

Further information can be found in the DiSH™ User's Guide (ENZ-71001-0001) on www.enzolifesciences.com.

REFERENCES

1. Price, P.W., McKinney, E.C., Wang, Y., Sasser, L.E., Kandasamy, M.K., Matsuuchi, L., Milcarek, C., Deal, R.B., Culver, D.G., and Meagher, R.B. (2009). Engineered cell surface expression of membrane immunoglobulin as a means to identify monoclonal antibody-secreting hybridomas. *J Immunol Methods* 343, 28-41.
2. Köhler G and Milstein C. Continuous culture of fused cells secreting antibody of predefined specificity. *Nature* 256:495-497, 1975
3. Kennett RH, McKearn TJ and Bechtol KB. Monoclonal antibodies. Plenum Press, New York 1980
4. Goding JW. Antibody production by hybridomas. *J. Immunol. Methods* 39:285-308, 1980
5. Sharon J, Morrison SL and Kabat EA. Detection of specific hybridoma clones by replica immunoadsorption of their secreted antibodies. *Proc. Natl. Acad.Sci.* 76:1420-1424, 1979
6. Pearson TW, Pinder M, Roelants G, Kar S, Lundin L, Mayor-Whitney KS and Hewett RS. Methods for derivation and detection of anti-parasite monoclonal antibodies. *J. Immunol. Methods.*34:141-154, 1980
7. Davis JM, Pennington JE, Kubler AM and Conscience JK. A simple, single-step technique for selecting and cloning hybridomas for the production of monoclonal antibodies. *J. Immunol. Methods* 50:161-171, 1982

For Research Use Only. This product and its use are the subject of US Patent Number 7,148,040.

Manufactured by 