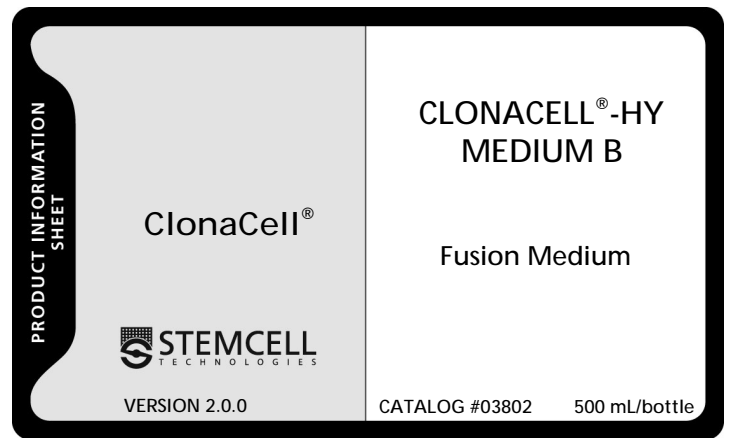


PRODUCT DESCRIPTION

This product contains gentamycin in Dulbecco's Modified Eagle's Medium (DMEM). This product is used to wash spleen cells and parental myeloma cells to remove all traces of serum prior to fusion of these cells.

STABILITY AND STORAGE

Product stable at -20°C until expiry date as indicated on label. Stable for four months when stored at 2 - 8°C. Product is sterility tested.



DIRECTIONS FOR USE

Thaw ClonaCell®-HY Medium B at 2 - 8°C overnight or at room temperature (15 - 25°C). Mix well. The product should not be thawed at 37°C or higher temperatures.

Refer to "ClonaCell®-HY Hybridoma Cloning Kit Technical Manual" for complete details available at www.stemcell.com/technical/manuals.aspx.

REFERENCES

1. Harlow E and Lane D. Monoclonal Antibodies, Chapter 6: Antibodies, a laboratory manual. Cold Spring Harbor Laboratory, Cold Sping Harbor, New York ISBN 0-87969-314-2, 1988
2. Köhler G and Milstein C. Continuous culture of fused cells secreting antibody of predefined specificity. Nature 256:495-497, 1975
3. Melchers F, Potter M and Warner NL. Lymphocyte hybridomas. Second workshop on "functional properties of tumours of T and B lymphocytes." Preface. Current Topics in Microbiology & Immunology 81:9-23, 1978
4. Kennett RH, McKearn TJ and Bechtol KB. Monoclonal antibodies. Plenum Press, New York 1980
5. Goding JW. Antibody production by hybridomas. J. Immunol. Methods 39:285-308, 1980
6. 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
7. 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
8. 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

