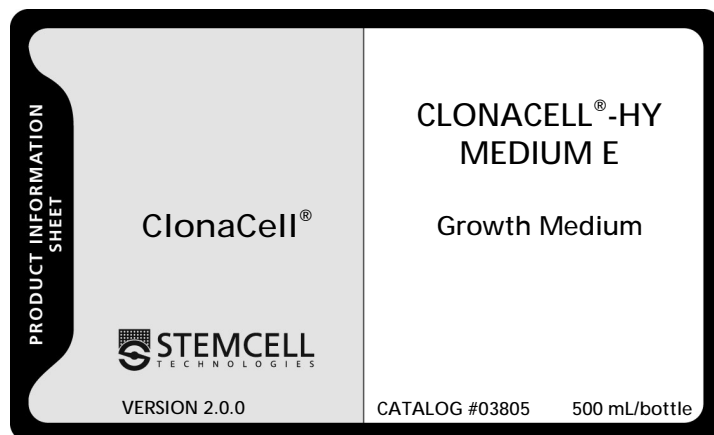


## PRODUCT DESCRIPTION

This product contains a mixture of pre-selected serum, Hypoxanthine Thymidine (HT), gentamycin and supplements in Dulbecco's Modified Eagle's Medium (DMEM) and has been formulated to promote the growth and expansion of hybridoma cell clones.

## STABILITY AND STORAGE

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



## DIRECTIONS FOR USE

Thaw ClonaCell®-HY Medium E 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](http://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

