

ES-Cult™ Basal Medium-A

Catalog # 05801

500 mL per bottle

This product has been pre-screened for use with other reagents of the ES-Cult™ product line for optimal performance for the *in vitro* pancreatic and neural differentiation of mouse embryonic stem (ES) cells.

PRODUCT DESCRIPTION:

This product contains: 1.55 g/L D-glucose, 0.5 mM L-glutamine, and 2.4 g/L Sodium Bicarbonate in Dulbecco's Modified Eagle's Medium (DME)/Ham's Nutrient Mixture F-12.

This product **does not** contain Antibiotics.

pH 7.2 –7.6

STABILITY / STORAGE:

Store at 2 -8°C.

Stable for one year from date of manufacture (refer to: Directions For Use).

Contents guaranteed sterile if seal is not tampered with.

DIRECTIONS FOR USE:

This product is recommended to be used in combination with ITS (Catalog #07151, #07155) or N2 Supplements (Catalog #07152, #07156) for *in vitro* pancreatic and neural differentiation of mouse ES cells. For complete instructions refer to the ES-Cult™ Technical Manuals on our website, www.stemcell.com.

This product contains the labile amino acid L-glutamine which has a half-life of approximately one month when stored at 2-8°C. For certain cell culture applications, medium stored for greater than two months following date of manufacture should be supplemented with additional L-glutamine. Add 1.25 mL of 200 mM L-glutamine (Catalog #07100) to 500 mL to achieve a final concentration of 0.5 mM.

If desired, add 5 mL of Penicillin and Streptomycin Solution (Catalog #07500) to each 500 mL bottle of this product to achieve a final concentration of 100 U/mL Penicillin and 100 µg/mL of Streptomycin. Following addition of antibiotics, medium should be used within one week.

**THIS REAGENT IS FOR LABORATORY USE ONLY.
IT IS NOT TO BE ADMINISTERED TO HUMANS.**