

ASCORBIC ACID

Catalog # 07157

100 milligrams per bottle

PRODUCT DESCRIPTION:

Ascorbic acid (C₆H₈O₆)
Vitamin C
Molecular weight 176.1 g/mol

STABILITY / STORAGE:

Store at room temperature. When dissolved in solution, protect from light and exposure to air to prevent oxidation.

DIRECTIONS FOR USE:

For ES-Cultä

This product has been developed to improve the yield of neural cells during the final stages of ES *in vitro* differentiation into neural cell types. It has been pre-screened for use with other reagents of the ES-Cult™ product line for optimal performance for the *in vitro* neural differentiation of murine ES cells.

Prepare a 0.02 M solution of ascorbic acid by adding 28.4 mL of sterile water to the amber bottle of ascorbic acid. Mix well, store at room temperature, protect from light and exposure to air. This provides a 100X stock solution and should be added freshly at each feed to the ES-Cult™ Basal Medium-A for a final concentration of 200 µM.

For more details the *in vitro* differentiation of murine ES cells to neural cell types, refer to protocols available on our website at www.stemcell.com.

For MesenCultä

This product is part of the Mesenchymal Stem Cell Osteogenic Stimulatory Supplement kit (catalog # 05404) used to stimulate the differentiation of human mesenchymal stem cells into osteogenic cells in culture.

This product should be added fresh each time to the Mesenchymal Stem Cell Osteogenic Stimulatory Supplements. A final concentration of 50 µg/mL is recommended.

For more details on the *in vitro* differentiation of human mesenchymal stem cells to osteogenic cells, refer to protocols available on our website at www.stemcell.com.

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

REFERENCES:

1. Lee SH, Lumelsky N, Studer L, Auerbach JM, and McKay RD. Efficient generation of midbrain neurons from mouse embryonic stem cells. *Nature Biotech.* 2000;**18**:675.