# **Cascade Biologics**™

invitrogen cell culture



# **SMDS**

## **Smooth Muscle Differentiation Supplement**

Cat. no. S-008-5

5 ml

### **Product Description**

Smooth Muscle Differentiation Supplement (SMDS) is a sterile, concentrated (100X) solution intended for use with basal Medium 231 for differentiation of normal human vascular smooth muscle cells. SMDS is an ionically-balanced supplement containing fetal bovine serum and heparin. Each 5 ml bottle of SMDS is the correct amount of differentiation supplement for a 500 ml bottle of basal Medium 231. When a 500 ml bottle of medium is supplemented with SMDS, the final concentrations of the components in the supplemented medium are: fetal bovine serum, 1% v/v; and heparin, 30  $\mu g/ml$ .

#### Intended Use

SMDS is intended for use in conjunction with basal Medium 231 to induce differentiation of normal human vascular smooth muscle cells. Vascular smooth muscle cells cultured in Medium 231 supplemented with Smooth Muscle Growth Supplement (SMGS, cat. # S-007-25) will proliferate. Upon transferring these proliferating cultures into Medium 231 supplemented with SMDS, the cells will undergo a cessation of growth, a change in cellular morphology from slender stellate cells to enlarged rectangular or triangular shaped cells, and will express increased amounts of smooth muscle  $\alpha\text{-actin}$  over the course of 2 to 8 days.

Reducing serum concentration and adding heparin to smooth muscle cell cultures are both commonly used procedures in differentiation studies of smooth muscle cells *in vitro*. For example see:

Newcomb, P.M. and Herman, I.M.; Pericyte growth and contractile phenotype: Modulation by endothelial-synthesized matrix and comparison with aortic smooth muscle. J. Cell Physiol 155: 385-393; 1993

Reilly, C.F., Fritze, L.M.S., Rosenburg, R.D.; Heparin inhibition of smooth muscle proliferation: a cellular site of action. J. Cell Physiol 129: 11-19; 1986

Castellot, J.J., Favreau, L.V., Karnovsky, M.J., Rosenberg, R.D.; Inhibition of vascular smooth muscle cell growth by endothelial cell-derived heparin. J. Biol Chem 257(19): 11256-11260; 1982.

This product is for research use only. Not for use in animals, humans, or diagnostic procedures.

Caution: If handled improperly, some components of this product may present a health hazard. Take appropriate precautions when handling this product, including the wearing of protective clothing and eyewear. Dispose of properly.

### Storage and Stability

SMDS is stored at -20°C at our facility and is shipped on dry ice. Upon receipt, the product should be stored at -20°C in a freezer that is <u>not</u> self-defrosting. When stored at -20°C the product is stable until the expiration date shown on the label.

### **Thawing**

To thaw, place the product in a 37°C water bath or overnight at 4°C. If thawed in a water bath, do not leave the product at 37°C after the product has thawed. For instructions on adding SMDS to basal Medium 231, please refer to the instructions that accompany the basal medium.

### For research use only.

## Cascade Biologics™

invitrogen cell culture



#### Limited Use Label License No. 5: Invitrogen Technology

The purchase of this product conveys to the buyer the non-trnsferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. Email: outlicensing@invitrogen.com.

©2009 Life Technologies Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.