

# Product Information



## Cell-Based Assay U-18666A

Item No. 10009869

### Contents

The vial contains 100 µl of 2.5 mM U-18666A in DMSO.

### Laboratory Procedures

Cell-Based Assay U-18666A is a cholesterol transport inhibitor and is provided for use as a positive control in the Cholesterol Cell-Based Detection Assay Kit, Cholesterol Uptake Cell-Based Assay Kit, and SREBP-2 Cell-Based Translocation Assay Kit (Item Nos. 10009779, 600440, and 10009239, respectively). The solution can be diluted up to 1:1,000 to treat experimental cells. This stock solution is good for one year when stored at -20°C.

### CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with both Cholesterol Cell-Based Detection Assay Kit, Cholesterol Uptake Cell-Based Assay Kit, and SREBP-2 Cell-Based Translocation Assay Kit from Cayman (Item Nos. 10009779, 600440, and 10009239, respectively). This product may not perform as described if used with other assay reagents or protocols.

### Related Products

Cholesterol Uptake Cell-Based Assay Kit - Item No. 600440 • SREBP-2 Cell-Based Translocation Assay Kit - Item No. 10009239 • Cholesterol Cell-Based Detection Assay Kit - Item No. 10009779

## Cayman Chemical

### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

### Phone

(800) 364-9897  
(734) 971-3335

### Fax

(734) 971-3640

### E-Mail

custserv@caymanchem.com

### Web

www.caymanchem.com

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/22/2011