

Product Information



Cell-Based Assay Soluble Epoxide Hydrolase Inhibitor (10 mM)

Item No. 600096

Contents: This vial contains 50 µl of 10 mM AUDA in DMSO

Stability: ≥1 year at -20°C

Laboratory Procedures

To make a Working Solution of Soluble Epoxide Hydrolase (sEH) Inhibitor, dilute 10 µl of this inhibitor with 500 µl of the diluted Assay Buffer. The diluted solution is stable for two hours on ice. If you are not using all of the sEH Inhibitor at one time, we recommend that you make small aliquots for long-term storage. The sEH Inhibitor, AUDA, is stable for one year at -20°C.

To test the Soluble Epoxide Hydrolase specificity of the samples, add 10 µl of this Working Solution along with 90 µl of sample supernatant to corresponding wells of a black plate. Mix with 100 µl of the diluted sEH Substrate Solution (Item No. 600095) and incubate at 37°C for 30 minutes.

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work with the sEH Cell-Based Assay Kit from Cayman (Item No. 600090). This product may not perform as described if used with other assay reagents or protocols.

Related Product

Soluble Epoxide Hydrolase Cell-Based Assay Kit - Item No. 600090

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, 01/26/2011