

Product Information



2-thio-PAF-AH Inhibitor Screening Assay (substrate)

Catalog No. 10004802

Laboratory Procedures

This vial contains an ethanolic solution of 2-thio PAF. 2-thio PAF, when stored as supplied, is stable for one year at -20°C. To prepare this substrate for use, evaporate the ethanolic solution to dryness under a gentle stream of nitrogen. Reconstitute the contents of the vial by vortexing with 12 ml of diluted PAF-AH Inhibitor Screening Assay Buffer (10X) (Catalog No. 10004800) to achieve a final concentration of 400 µM. Make sure to vortex until the substrate solution becomes clear. The reconstituted substrate is stable for two weeks at -20°C. Two substrate vials are required to perform the entire PAF Acetylhydrolase Inhibitor Screening Assay Kit (Catalog No. 10004380).

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 the PAF Acetylhydrolase Inhibitor Screening Assay Kit from Cayman Chemical (Catalog No. 10004380). This product may not perform as described if used with other assay reagents or protocols.

Related Products

PAF Acetylhydrolase Inhibitor Screening Assay Kit - Cat. No. 10004380 • PAF-AH Inhibitor Screening Assay Buffer (10X) - Cat. No. 10004800 • PAF-AH Inhibitor Screening Assay DTNB - Cat. No. 10004801 • Human Plasma PAF-AH Inhibitor Screening Assay - Cat. No. 10004803 • 96-WellCover Sheets - Cat. No. 400012 • 96-Well Plate (Colorimetric Assay) - Cat. No. 400014

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, 11/20/2009