

Product Information



cPLA₂ Assay Buffer

Item No. 765030

Contents

The vial contains 30 ml of 160 mM Hepes buffer (pH 7.4) containing 20 mM CaCl₂, 300 mM sodium chloride, 8 mM Triton X-100, 60% glycerol, and 2 mg/ml BSA. Store at -20°C. When stored under these conditions, the concentrated buffer is stable for at least one year. This concentrated buffer is used directly (without dilution) for the reconstitution of Arachidonoyl Thio-PC (substrate) (Item No. 765034).

Laboratory Procedures

The Assay Buffer is diluted before assaying for calcium dependent cPLA₂ activity, using the following procedure. Dilute 10 ml of the Assay Buffer concentrate with 10 ml of HPLC-grade (or equivalent) water. This final Assay Buffer contains 80 mM Hepes (pH 7.4) containing 10 mM CaCl₂, 150 mM sodium chloride, 4 mM Triton X-100, 30% glycerol, and 1 mg/ml BSA and should be used for the dilution of samples prior to assaying for cPLA₂ activity. This diluted Assay Buffer is stable for at least one month if stored at 4°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 the cPLA₂ Assay Kit from Cayman (Item No. 765021). This product may not perform as described if used with other assay reagents or protocols.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/765030

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, 01/31/2013

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