

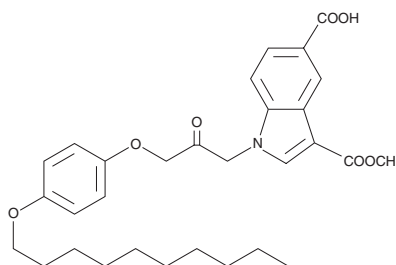
Product Information



CAY10502

Item No. 10008657

Formal Name: 3-methylhydrogen-1-[3-(4-decyloxy phenoxy)-2-oxopropyl]indole-3,5-dicarboxylate
MF: C₃₀H₃₇NO₇
FW: 523.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10502 be stored as supplied at -20°C. It will be stable for at least two years.

CAY10502 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10502 in an organic solvent purged with an inert gas. CAY10502 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CAY10502 in DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

CAY10502 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10502 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10502 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Phospholipase A₂ (PLA₂) catalyzes the hydrolysis of phospholipids at the *sn*-2 position leading to the production of lysophospholipids and free fatty acids. Calcium-dependent cytosolic PLA₂α (cPLA₂α) is a 85 kDa enzyme that plays a key role in the arachidonic cascade and the inflammatory response associated with this metabolic pathway.¹ CAY10502 is a potent inhibitor of cPLA₂α with an IC₅₀ value of 4.3 nM for the purified enzyme from human platelets.² It inhibits arachidonic acid mobilization from A23187-stimulated or TPA-stimulated human platelets with IC₅₀ values of 570 and 0.9 nM, respectively.²

References

- Schaloske, R.H. and Dennis, E.A. The phospholipase A₂ superfamily and its group numbering system. *Biochem. Biophys. Acta* **1761**, 1246-1259 (2006).
- Ludwig, J., Bovens, S., Brauch, C., *et al.* Design and synthesis of 1-indol-1-yl-propan-2-ones as inhibitors of human cytosolic phospholipase A₂α. *J. Med. Chem.* **49**, 2611-2620 (2006).

Related Products

RSC-3388 - Item No. 10007784 • CAY10641 - Item No. 10662

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, 03/07/2011