

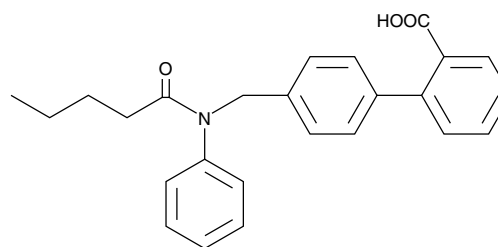
Product Information



CAY10583

Item No. 10012424

CAS Registry No.: 862891-27-8
Formal Name: 4'-[[[(1-oxopentyl)phenylamino]methyl]-[1,1'-biphenyl]-2-carboxylic acid
MF: C₂₅H₂₅NO₃
FW: 387.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10583 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10583 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10583 in an organic solvent purged with an inert gas. CAY10583 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10583 in ethanol is approximately 0.25 mg/ml and approximately 10 mg/ml in DMSO and DMF.

CAY10583 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10583 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10583 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Leukotriene B₄ (LTB₄) promotes a number of leukocyte functions including aggregation, stimulation of ion fluxes, superoxide anion production, chemotaxis, and chemokinesis. Two G protein-coupled receptors, BLT₁ and BLT₂ mediate the actions of LTB₄. CAY10583 is a potent, selective full agonist for BLT₂ with an EC₅₀ value of 20 nM.¹ It does not displace [³H]-LTB₄ binding to BLT₁ at concentrations >1 μM and activates BLT₂ at significantly lower concentrations than LTB₄ (EC₅₀ = 170 nM).¹ CAY10583 dose dependently increases intracellular calcium stores and induces ERK phosphorylation in Chinese hamster ovary (CHO) cells expressing BLT₂ with no effect on CHO cells expressing BLT₁.¹

Reference

1. Iizuka, Y., Yokomizo, T., Terawaki, K., *et al.* Characterization of a mouse second leukotriene B₄ receptor, mBLT₂. BLT₂-dependent ERK activation and cell migration of primary mouse keratinocytes. *J. Biol. Chem.* **280(26)**, 24816-24823 (2005).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10012424

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

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 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, 05/02/2014

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