

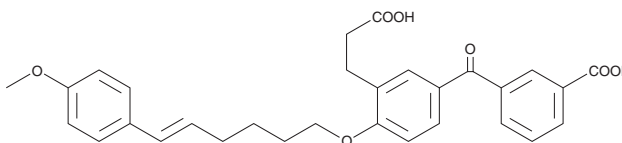
Product Information



LY223982

Catalog No. 10010024

CAS Registry No.: 117423-74-2
Formal Name: 5-(3-carboxybenzoyl)-2-
[[[(5E)-6-(4-methoxyphenyl)-
5-hexenyl]oxy]-
benzenepropanoic acid
Synonyms: CGS 23131, SKF 107324
MF: C₃₀H₃₀O₇
FW: 502.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that LY223982 be stored as supplied at -20°C. It should be stable for at least two years.

LY223982 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY223982 in an organic solvent purged with an inert gas. LY223982 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LY223982 in ethanol is approximately 0.1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

LY223982 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY223982 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LY223982 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.

Leukotriene B₄ (LTB₄) is a dihydroxy fatty acid derived from the 5-lipoxygenase (5-LO) pathway of arachidonic acid metabolism and is an important mediator of the inflammatory process. The benzophenone dicarboxylic acid derivative LY223982 is a potent BLT₁ receptor antagonist. It inhibits the specific binding of radiolabeled-LTB₄ to isolated human neutrophils with an IC₅₀ value of 13.2 nM.¹ LY223982 inhibits the LTB₄-induced aggregation of guinea pig and human neutrophils with IC₅₀ values of 74 and 100 nM, respectively.² However, concentrations of LY223982 up to 10 μM do not inhibit binding of LTB₄ to human BLT₁ or BLT₂ expressed in CHO cells.³

References

1. Gapinski, D.M., Mallett, B.E., Froelich, L.L., *et al.* Benzophenone dicarboxylic acid antagonists of leukotriene B₄. 2. Structure-activity relationships of the lipophilic side chain. *J. Med. Chem.* **33**, 2807-2813 (1990).
2. Jackson, W.T., Boyd, R.J., Froelich, L.L., *et al.* Specific inhibition of leukotriene B₄-induced neutrophil activation by LY223982. *J. Pharmacol. Exp. Ther.* **263**(3), 1009-1014 (1992).
3. Yokomizo, T., Kato, K., Hagiya, H., *et al.* Hydroxyeicosanoids bind to and activate the low affinity leukotriene B₄ receptor, BLT₂. *J. Biol. Chem.* **276**(15), 12454-12459 (2001).

Related Product

LY293111 - Cat. No. 10009768

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, 11/01/2010

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