

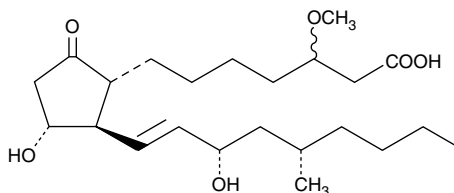
Product Information



3-methoxy Limaprost

Item No. 10008445

CAS Registry No.: 1224443-96-2
Formal Name: 3-methoxy-11 α ,15S-dihydroxy-17S,20-dimethyl-9-oxo-prosta-13E-en-1-oic acid
Synonyms: 3-methoxy-2,3-dihydro-Limaprost, 3-methoxy-17 α ,20-dimethyl-Prostaglandin E₁
MF: C₂₃H₄₀O₆
FW: 412.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 3-methoxy limaprost be stored as supplied at -20°C. It should be stable for at least one year.

3-methoxy Limaprost is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 3-methoxy limaprost in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 3-methoxy limaprost is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 3-methoxy limaprost in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Limaprost is an analog of prostaglandin E₁ (PGE₁) with structural modifications intended to give it a prolonged half-life and greater potency. Limaprost is an inhibitor of ADP and collagen-induced platelet aggregation and is 10-1,000 times more potent than PGE₁ as an inhibitor of platelet adhesiveness measured *in vitro*.¹ 3-methoxy Limaprost results from the Michael addition of methanol to the limaprost 2,3 double bond. The resulting novel compound is likely to retain FP receptor agonist activity; however there are no published studies of the pharmacological properties of 3-methoxy limaprost.

Reference

1. Tsuboi, T., Hatano, N., Nakatsuji, K., *et al.* Pharmacological evaluation of OP 1206, a prostaglandin E₁ derivative, as an antianginal agent. *Arch. Int. Pharmacodyn.* **247**, 89-102 (1980).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008445

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, 08/05/2013

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