

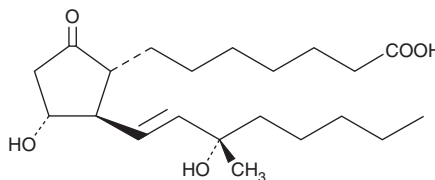
PRODUCT INFORMATION



15(S)-15-methyl Prostaglandin E₁

Item No. 13730

CAS Registry No: 35700-26-6
Formal Name: 9-oxo-11 α ,15S-dihydroxy-15-methyl-prost-13E-en-1-oic acid
Synonym: 15(S)-15-methyl PGE₁,
MF: C₂₁H₃₆O₅
FW: 368.5
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 15(S)-15-methyl prostaglandin E₁ (15(S)-15-methyl PGE₁) be stored as supplied at -20°C. It should be stable for at least two years.

15(S)-15-methyl PGE₁ 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 15(S)-15-methyl PGE₁ in these solvents is approximately 50 mg/ml. 15(S)-15-methyl PGE₁ is stable for at least six months in these solvents if stored at -20°C.

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 15(S)-15-methyl PGE₁ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 15(S)-15-methyl PGE₁ in PBS (pH 7.2) is approximately 1.6 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15(S)-15-methyl PGE₁ is a metabolically stable synthetic analog of PGE₁ (Item No. 13010). In sharp contrast to the bronchodilatory effects of PGE₁, 15(S)-15-methyl PGE₁ is a weak constrictor of human respiratory tract smooth muscle.¹

Reference

1. Karim, S.M.M., Adaikan, P.G., and Kottegoda, S.R. Prostaglandins and human respiratory tract smooth muscle: Structure activity relationship. *Adv. Prostaglandin Thromboxane Res.* **7**, 969-980 (1980).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/11/2015

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM