

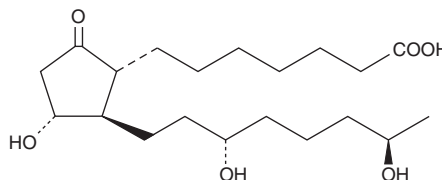
PRODUCT INFORMATION



13,14-dihydro-19(R)-Prostaglandin E₁

Item No. 13660

Formal Name: 9-oxo-11 α ,15S,19R-trihydroxy-prostan-1-oic acid
Synonyms: 13,14-dihydro-19(R)-PGE₁, 19(R)-hydroxy PGE₀
MF: C₂₀H₃₆O₆
FW: 372.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 13,14-dihydro-19(R)-hydroxy prostaglandin E₁ (13,14-dihydro-19(R)-hydroxy PGE₁) be stored as supplied at -20°C. It should be stable for at least one year.

13,14-dihydro-19(R)-hydroxy PGE₁ is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 13,14-dihydro-19(R)-hydroxy PGE₁ in these solvents is approximately 50 mg/ml. 13,14-dihydro-19(R)-hydroxy 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 13,14-dihydro-19(R)-hydroxy PGE₁ is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 13,14-dihydro-19(R)-hydroxy 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

13,14-dihydro-19(R)-hydroxy PGE₁ is the theoretical metabolite of 13,14-dihydro PGE₁ (Item No. 13610) by ω -1 hydroxylase. There are no known reports on the biosynthesis or biological activity of 13,14-dihydro-19(R)-hydroxy PGE₁.

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