

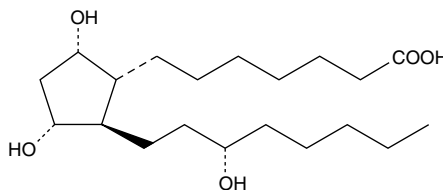
Product Information



13,14-dihydro Prostaglandin F_{1α}

Item No. 15610

CAS Registry No.: 20592-20-5
Formal Name: 9α,11α,15S-trihydroxy-prostan-1-oic acid
Synonym: Prostaglandin F_α
MF: C₂₀H₃₈O₅
FW: 358.5
Purity: ≥99%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

13,14-dihydro Prostaglandin F_{1α} (13,14-dihydro PGF_{1α}) is a metabolite of PGF_{1α}. For long term storage, we suggest that 13,14-dihydro PGF_{1α} be stored as supplied at -20°C. It will be stable for at least two years.

13,14-dihydro PGF_{1α} is supplied as a crystalline solid. 13,14-dihydro PGF_{1α} is soluble in organic solvents such as ethanol, DMSO, or dimethyl formamide. The solubility of 13,14-dihydro PGF_{1α} in these solvents is approximately 50 mg/ml. 13,14-dihydro PGF_{1α} 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. Organic solvent-free aqueous solutions of 13,14-dihydro PGF_{1α} can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 13,14-dihydro PGF_{1α} in PBS (pH 7.2) is approximately 2.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

13,14-dihydro PGF_{1α}, 13,14-dihydro-15-keto PGF_{1α}, and 13,14-dihydro-15-keto PGF_{2α} are initial metabolites of PGF_{1α} and PGF_{2α}.¹ The measurement of these metabolites in cell culture is used to assess total PGF metabolism.¹

Reference

1. Ritz, E.M., Boto, W.O., Stylos, W.A. Measurement of initial prostaglandin F metabolites in medium of BALB/c 3T3 and SV3T3 mouse fibroblast cultures. *Biochem. Biophys. Res. Commun.* **63**, 179-186 (1975).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/15610

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, 05/31/2013

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