

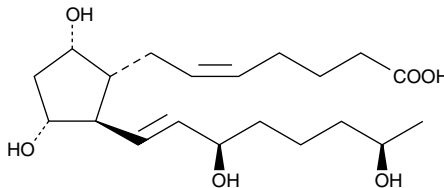
Product Information



15(R),19(R)-hydroxy Prostaglandin F_{2α}

Item No. 16920

CAS Registry No.: 1224444-23-8
Formal Name: 9α,11α,15R,19R-tetrahydroxy-prosta-5Z,13E-dien-1-oic acid
Synonym: 15(R),19(R)-hydroxy PGF_{2α}
MF: C₂₀H₃₄O₆
FW: 370.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 15(R),19(R)-hydroxy prostaglandin F_{2α} (15(R),19(R)-hydroxy PGF_{2α}) be stored as supplied at -20°C. It should be stable for at least one year.

15(R),19(R)-hydroxy PGF_{2α} 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 15(R),19(R)-hydroxy PGF_{2α} in these solvents is approximately 100 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 15(R),19(R)-hydroxy PGF_{2α} is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15(R),19(R)-hydroxy PGF_{2α} in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

19(R)-Hydroxylated PGs occur in μg/ml concentrations in the semen of certain mammalian species, especially primates. In the case of humans, the compounds are primarily of the PGE series, and the hydroxyl stereochemistry is 15(S),19(R).¹ 19(R)-Hydroxylated PGs are also found in the seminal plasma of marsupials, where F-type compounds of the 1 and 2-series predominate.² The 15(R)-hydroxy epimer of these 19-hydroxylated PGs is the inverse or 'unnatural' isomer at C-15. The biological role of 19(R)-hydroxylated PGs is not well established. In the F-series, 19(R)-hydroxylation is associated with a significant loss of receptor-mediated biological activity in some assays.³

References

1. Cooper, I. and Kelly, R.W. The measurement of E and 19-hydroxy E prostaglandins in human seminal plasma. *Prostaglandins* **10**, 507-514 (1975).
2. Marley, P.B., Rodger, J.C., White, I.G., *et al.* 19-Hydroxylated prostaglandins in the semen of the marsupial *Trichosurus vulpecula* (brush-tailed possum). *Comp. Biochem. Physiol.* **70B**, 619-621 (1981).
3. Woodward, D.F., Protzman, C.E., Krauss, A.H.P., *et al.* Identification of 19(R)-OH prostaglandin E₂ as a selective prostanoid EP2-receptor agonist. *Prostaglandins* **46**, 371-383 (1993).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16920

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, 02/06/2014

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