

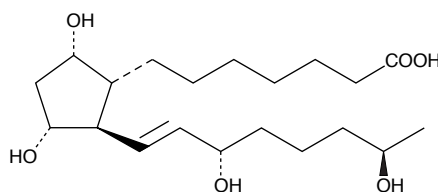
# Product Information



## 19(R)-hydroxy Prostaglandin F<sub>1α</sub>

Catalog No. 15910

**CAS Registry No:** 81371-59-7  
**Formal Name:** 9α,11α,15S,19R-tetrahydroxy-prost-13E-en-1-oic acid  
**MF:** C<sub>20</sub>H<sub>36</sub>O<sub>6</sub>  
**FW:** 372.5  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A solution in ethanol



### Laboratory Procedures

For long term storage, we suggest that 19(R)-hydroxy Prostaglandin F<sub>1α</sub> (19(R)-hydroxy PGF<sub>1α</sub>) be stored as supplied at -20°C. It will be stable for at least two years.

19(R)-hydroxy PGF<sub>1α</sub> 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 19(R)-hydroxy PGF<sub>1α</sub> in these solvents is approximately 50 mg/ml. 19(R)-hydroxy PGF<sub>1α</sub> 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 19(R)-hydroxy PGF<sub>1α</sub> is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 19(R)-hydroxy PGF<sub>1α</sub> in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

19(R)-hydroxy PGF<sub>1α</sub> is an ω-1 hydroxylase metabolite of PGF<sub>1α</sub> that has been identified in the semen of humans and marsupials.<sup>1,2</sup> There are no published reports on the biological activity of 19(R)-hydroxy PGF<sub>1α</sub>.

### References

1. Taylor, P.L. and Kelly, R.W. The occurrence of 19-hydroxy F prostaglandins in human semen. *FEBS Lett.* **57**, 22-25 (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).

### Related Product

19(R)-hydroxy Prostaglandin E<sub>1</sub> - Cat. No. 13910 • 19(R)-hydroxy Prostaglandin E<sub>2</sub> - Cat. No. 14910 • Prostaglandin F<sub>1α</sub> - Cat. No. 15010 • 15(R),19(R)-hydroxy Prostaglandin F<sub>1α</sub> - Cat. No. 15910 • 19(R)-hydroxy Prostaglandin F<sub>2α</sub> - Cat. No. 16910

**WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### MATERIAL 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 Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at 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, 04/04/2006

### 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