

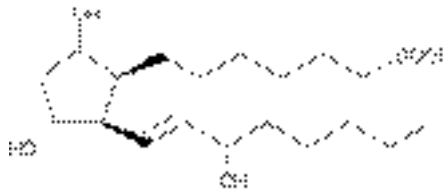
# Product Information



## 8-*iso* Prostaglandin F<sub>1α</sub>

Item No. 15350

**CAS Registry No.:** 26771-96-0  
**Formal Name:** 9α,11α,15S-trihydroxy-(8β)-prost-13E-en-1-oic acid  
**Synonym:** 8-*epi* Prostaglandin F<sub>1α</sub>  
**MF:** C<sub>20</sub>H<sub>36</sub>O<sub>5</sub>  
**FW:** 356.5  
**Purity:** ≥99%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

8-*iso* Prostaglandin F<sub>1α</sub> (8-*iso* PGF<sub>1α</sub>) is an isoprostane that was first identified in human semen.<sup>1</sup> For long term storage, we suggest that 8-*iso* PGF<sub>1α</sub> be stored as supplied at -20°C. It will be stable for at least two years.

8-*iso* PGF<sub>1α</sub> is supplied as a crystalline solid. A stock solution may be made by dissolving the 8-*iso* PGF<sub>1α</sub> in an organic solvent. 8-*iso* PGF<sub>1α</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 8-*iso* PGF<sub>1α</sub> in these solvents is approximately 50 mg/ml. 8-*iso* PGF<sub>1α</sub> will be 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 8-*iso* PGF<sub>1α</sub> can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 8-*iso* PGF<sub>1α</sub> in PBS (pH 7.2) is approximately 2.0 mg/ml. We do not recommend storing the aqueous solution for more than one day.

8-*iso* PGF<sub>1α</sub> is a member of the isoprostane family, which are eicosanoids of non-cyclooxygenase origin. 8-*iso* PGF<sub>1α</sub> is found in human semen along with its 19-hydroxy congener at 5-10 µg/ml of seminal plasma.<sup>2</sup>

### References

1. Svanborg, K., Bygdeman, M., and Eneroth, P. The F and 19-hydroxy F prostaglandins and their 8-isomers in human seminal plasma: Data on chromatography and mass spectrometry. *Biomedical Mass Spectrometry* **10**, 495-498 (1983).
2. Taylor, P.L. The 8-isoprostaglandins: Evidence for eight compounds in human semen. *Prostaglandins* **17**, 259-267 (1979).

### Related Product

19(R)-hydroxy Prostaglandin F<sub>1α</sub> - Item No. 15910

**WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.**

#### 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, 10/03/2011

### 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