

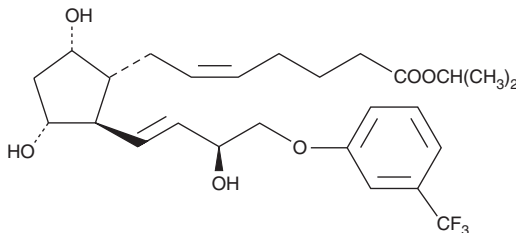
# Product Information



## 15(S)-Fluprostenol isopropyl ester

Item No. 16788

**CAS Registry No.:** 1420791-14-5  
**Formal Name:** 9 $\alpha$ ,11 $\alpha$ ,15S-trihydroxy-16-(3-trifluoromethyl)phenoxy)-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-oic acid, isopropyl ester  
**Synonym:** 15(S)-Flu-Ipr  
**MF:** C<sub>26</sub>H<sub>35</sub>F<sub>3</sub>O<sub>6</sub>  
**FW:** 500.6  
**Purity:**  $\geq$ 98%  
**Stability:**  $\geq$ 2 years at -20°C  
**Supplied as:** A solution in ethanol  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 222, 280 nm



### Laboratory Procedures

For long term storage, we suggest that 15(S)-fluprostenol isopropyl ester (15(S)-Flu-Ipr) be stored as supplied at -20°C. It should be stable for at least two years.

15(S)-Flu-Ipr 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 15(S)-Flu-Ipr in these solvents is approximately 12 mg/ml. 15(S)-Flu-Ipr is stable for at least six months in these solvents if stored at -20°C.

15(S)-Flu-Ipr is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 15(S)-Flu-Ipr should be diluted with the aqueous buffer of choice. 15(S)-Flu-Ipr has a solubility of 1 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

15(S)-Flu-Ipr is the unnatural C-15 epimer of Travoprost. Travoprost is the Alcon trade name for Flu-Ipr an F-series prostaglandin analog which has been approved for use as an ocular hypotensive drug.<sup>1</sup> Fluprostenol isopropyl ester is a prodrug which is converted by esterase enzymatic activity in the cornea to yield the corresponding free acid. The free acid, fluprostenol, is a potent FP receptor agonist.<sup>2</sup> In human and animal models of glaucoma, FP receptor agonist activity corresponds very closely with intraocular hypotensive activity. Although data is currently not available for the activity of 15(S)-Flu-Ipr, inversion of the stereochemistry at the 15 position of ocular hypotensive prostaglandins generally lowers the potency approximately 100-fold.<sup>3</sup>

### References

1. Sorbera, L.A. and Castañer, J. Travoprost. *Drugs of the Future* **25**, 41-45 (2000).
2. Abramovitz, M., Adam, M., Boie, Y., *et al.* The utilization of recombinant prostanoid receptors to determine the affinities and selectivities of prostaglandins and related analogs. *Biochim. Biophys. Acta* **1483**, 285-293 (2000).
3. Stjerschantz, J. and Resul, B. Phenyl substituted prostaglandin analogs for glaucoma treatment. *Drugs of the Future* **17**, 691-704 (1992).

**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, 07/02/2015

### 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