

PRODUCT INFORMATION

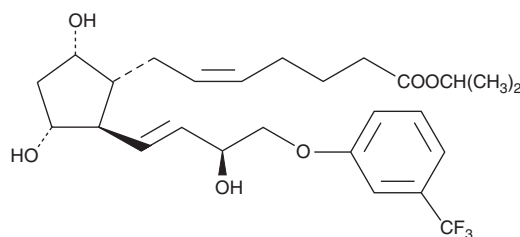


15(S)-Fluprostenol isopropyl ester

Item No. 16788

CAS Registry No.: 1420791-14-5
Formal Name: (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(1E,3S)-3-hydroxy-4-[3-(trifluoromethyl)phenoxy]-1-buten-1-yl]cyclopentyl]-5-heptenoic acid, 1-methylethyl ester

Synonym: 15(S)-Flu-lpr
MF: C₂₆H₃₅F₃O₆
FW: 500.6
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 280 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

15(S)-Fluprostenol isopropyl ester 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)-fluprostenol isopropyl ester in these solvents is approximately 30, 12, and 15 mg/ml, respectively.

15(S)-Fluprostenol isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 15(S)-fluprostenol isopropyl ester should be diluted with the aqueous buffer of choice. 15(S)-Fluprostenol isopropyl ester 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)-Fluprostenol isopropyl ester is an isomer of the prostaglandin F_{2α} analog fluprostenol isopropyl ester (Item No. 16769).¹ It is a potential prodrug to 15(S)-fluprostenol (Item No. 16787), which could act as an agonist at FP receptors albeit at a lower potency than 15(R) epimer and FP receptor agonist fluprostenol (Item No. 16768). 15(S)-Fluprostenol isopropyl ester is a potential impurity found in commercial preparations of fluprostenol isopropyl ester.²

References

1. Sorbera, L.A. and Castañer, J. Travoprost. *Drugs Future* **25**(1), 41-45 (2000).
2. Dams, I., Chodyński, M., Krupa, M., *et al.* A novel convergent synthesis of the potent antiglaucoma agent travoprost. *Tetrahedron* **69**(5), 1634-1648 (2013).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/11/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM