

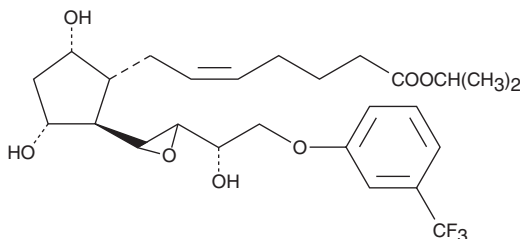
PRODUCT INFORMATION



13,14-epoxy Fluprostenol isopropyl ester

Item No. 13679

CAS Registry No.: 2557327-99-6
Formal Name: 7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(2R)-3-[(1S)-1-hydroxy-2-[3-(trifluoromethyl)phenoxy]ethyl]-2-oxiranyl]cyclopentyl]-5-heptenoic acid, 1-methylethyl ester
Synonym: 13,14-epoxy Travoprost
MF: C₂₆H₃₅F₃O₇
FW: 516.6
Purity: ≥95% (epoxide mixture)
UV/Vis.: λ_{max}: 222, 276 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

13,14-epoxy Fluprostenol isopropyl ester is supplied as methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 13,14-epoxy fluprostenol isopropyl ester in DMSO is approximately 5 mg/ml and approximately 30 mg/ml in ethanol and DMF.

13,14-epoxy Fluprostenol isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 13,14-epoxy fluprostenol isopropyl ester should be diluted with the aqueous buffer of choice. 13,14-epoxy Fluprostenol isopropyl ester has a solubility of 0.50 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

Fluprostenol isopropyl ester (Item No. 16769) is a potent F-series prostaglandin receptor agonist that finds clinical use as an ocular hypotensive agent for the treatment of glaucoma.^{1,2} 13,14-epoxy Fluprostenol isopropyl ester is an impurity generated in the production of fluprostenol isopropyl ester. This product is a 4:96 diastereomeric mixture of epoxides.

References

1. Sorbera, L.A. and Castañer, J. Travoprost. *Drugs of the Future* **25**, 41-45 (2000).
2. Hellberg, M.R., Sallee, V.L., Mclaughlin, M.A., et al. Preclinical efficacy of travoprost, a potent and selective FP prostaglandin receptor agonist. *J. Ocul. Pharmacol. Ther.* **17(5)**, 421-432 (2001).

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, 03/13/2024

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