

PRODUCT INFORMATION



Fluprostenol isopropyl ester-d₄

Item No. 9000936

Formal Name: 7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[[[(1E,3R)-3-hydroxy-4-[3-(trifluoromethyl)phenoxy]-1-buten-1-yl]cyclopentyl]-5Z-heptenoic-d₄ acid, 1-methylethyl ester

Synonyms: Flu-lpr-d₄, 16-m-trifluoromethylphenoxy tetranor PGF_{2α} isopropyl ester-d₄, Travoprost-d₄

MF: C₂₆H₃₁D₄F₃O₆

FW: 504.6

Chemical Purity: ≥98% (Fluprostenol isopropyl ester)

Deuterium

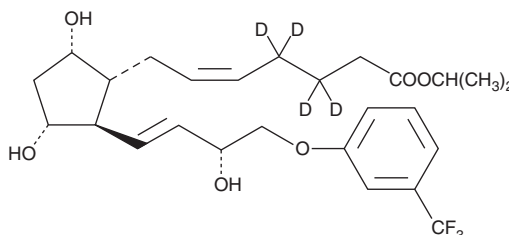
Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

UV/Vis.: λ_{max}: 222, 276 nm

Supplied as: A 500 µg/ml 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

Fluprostenol isopropyl ester-d₄ is intended for use as an internal standard for the quantification of fluprostenol isopropyl ester (Item No. 16769) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Fluprostenol isopropyl ester-d₄ is supplied as a solution in 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 purged with an inert gas can be used. The solubility of fluprostenol isopropyl ester-d₄ in these solvents is approximately 25, 20, and 30 mg/ml, respectively.

Description

Fluprostenol isopropyl ester is an analog of prostaglandin F_{2α} (PGF_{2α}; Item Nos. 16010 | 16020) and an isopropyl ester prodrug form of (+)-fluprostenol (Item No. 16768).^{1,2} Fluprostenol isopropyl ester is an FP receptor agonist, inducing phosphoinositide turnover in HEK293 cells expressing the human ocular FP receptor with an EC₅₀ value of 40.2 nM.³ Topical application of fluprostenol isopropyl ester (0.01, 0.03, and 0.1 µg) induces miosis in conscious cats in a dose-dependent manner.² It reduces intraocular pressure in a cynomolgus monkey model of ocular hypertension when administered topically at doses of 0.1 and 0.3 µg twice per day. Formulations containing fluprostenol isopropyl ester have been used in the treatment of open-angle glaucoma and ocular hypertension.

References

1. Sorbera, L.A. and Castañer, J. Travoprost. *Drugs Future* **25**(1), 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).
3. Sharif, N.A., Kelly, C.R., Crider, J.Y., et al. Ocular hypotensive FP prostaglandin (PG) analogs: PG receptor subtype binding affinities and selectivities, and agonist potencies at FP and other PG receptors in cultured cells. *J. Ocul. Pharmacol. Ther.* **19**(6), 501-515 (2003).

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, 11/24/2025

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