

# PRODUCT INFORMATION



## 13(R),14(R)-epoxy Fluprostenol isopropyl ester

Item No. 11169

**CAS Registry No.:** 2557329-36-7  
**Formal Name:** 7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(2R,3R)-3-[(1S)-1-hydroxy-2-[3-(trifluoromethyl)phenoxy]ethyl]-2-oxiranyl]cyclopentyl]-(5Z)-5-heptenoic acid, 1-methylethyl ester

**Synonyms:** Epoxy Derivative 1,  
13(R),14(R)-epoxy Travoprost

**MF:** C<sub>26</sub>H<sub>35</sub>F<sub>3</sub>O<sub>7</sub>

**FW:** 516.6

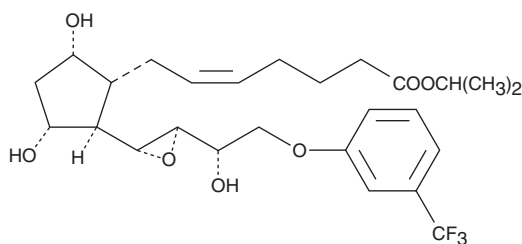
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 222, 276 nm

**Supplied as:** A solution in methyl acetate

**Storage:** -20°C

**Stability:** ≥1 year



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

### Laboratory Procedures

13(R),14(R)-epoxy Fluprostenol isopropyl ester 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 (DMF) purged with an inert gas can be used. The solubility of 13(R),14(R)-epoxy Fluprostenol isopropyl ester in ethanol and DMF is approximately 30 mg/ml and approximately 5 mg/ml in DMSO.

13(R),14(R)-epoxy Fluprostenol isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 13(R),14(R)-epoxy fluprostenol isopropyl ester should be diluted with the aqueous buffer of choice. 13(R),14(R)-epoxy Fluprostenol isopropyl ester has a solubility of 0.5 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

13,14-epoxy Fluprostenol isopropyl ester (Item No. 13679) is an impurity generated in the production of fluprostenol isopropyl ester. This product is a chiral enantiomer of the epoxide. The pharmacology of 13(R),14(R)-epoxy fluprostenol isopropyl ester has not been studied extensively to date.

#### 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, 05/20/2022

#### 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