

# PRODUCT INFORMATION

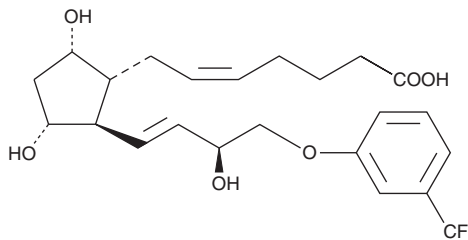


## 15(S)-Fluprostenol

Item No. 16787

**CAS Registry No.:** 54276-24-3  
**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

**MF:** C<sub>23</sub>H<sub>29</sub>F<sub>3</sub>O<sub>6</sub>  
**FW:** 458.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 277 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 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 15(S)-fluprostenol in these solvents is approximately 20 and 30 mg/ml, respectively.

15(S)-Fluprostenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 15(S)-fluprostenol should be diluted with the aqueous buffer of choice. The solubility of 15(S)-fluprostenol in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

15(S)-Fluprostenol is an isomer of the FP receptor agonist fluprostenol (Item No. 16768).<sup>1</sup> It is a potential active metabolite of 15(S)-fluprostenol isopropyl ester (Item No. 16788), which could act as an agonist at FP receptors albeit at a lower potency than the 15(R) epimer fluprostenol.

### Reference

1. Niyaz, M., Khan, M.S., and Mudassar, S. Hedgehog signaling: An achilles' heel in cancer. *Transl. Oncol.* **12(10)**, 1334-1344 (2019).

#### 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, 12/13/2023

#### 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