

PRODUCT INFORMATION



5-*trans* Fluprostenol

Item No. 16776

CAS Registry No.: 57968-83-9
Formal Name: [1R-[1 α (E),2 β (1E,3R*),3 α ,5 α]-7-(3,5-dihydroxy-2-[3-hydroxy-4-[3-(trifluoromethyl)phenoxy]-1-butenyl]cyclopentyl]-5-heptenoic acid

Synonym: 5,6-*trans* Fluprostenol

MF: C₂₃H₂₉F₃O₆

FW: 458.5

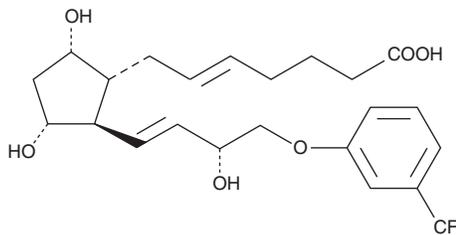
Purity: \geq 95%

UV/Vis.: λ_{max} : 222, 277 nm

Supplied as: A solution in ethanol

Storage: -20°C

Stability: \geq 2 years



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

Laboratory Procedures

5-*trans* 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 5-*trans* fluprostenol in these solvents is approximately 100 mg/ml.

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

Description

5-*trans* Fluprostenol is a potential impurity in commercial preparations of the FP receptor agonist fluprostenol (Item No. 16768).¹

Reference

1. Vajda, E., Hortobágyi, I., Lászlófi, I., et al. New process for the preparation of high purity prostaglandins. *Chinoin Gyógyszer és Vegyészeti Termékek Gyára Zrt.* **WO2015136317A1** (2015).

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, 11/30/2023

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