

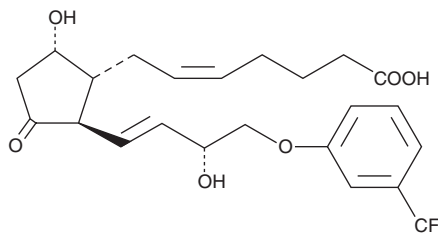
PRODUCT INFORMATION



11-keto Fluprostenol

Item No. 16783

CAS Registry No.: 62145-07-7
Formal Name: 7-[(1R,2R,5S)-5Z-hydroxy-2-[(1E,3R)-3-hydroxy-4-[3-(trifluoromethyl)phenoxy]-1-buten-1-yl]-3-oxocyclopentyl]-5-heptenoic acid
Synonym: Fluprostenol Prostaglandin D₂
MF: C₂₃H₂₇F₃O₆
FW: 456.5
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 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

11-keto Fluprostenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 11-keto fluprostenol should be diluted with the aqueous buffer of choice. The solubility of 11-keto fluprostenol in PBS (pH 7.2) is approximately 29 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

11-keto Fluprostenol is an analog of prostaglandin D₂ (PGD₂) with structural modifications intended to give it a prolonged half-life and greater potency. Fluprostenol is a well-studied, potent analog of PGF_{2α} and acts primarily through the FP receptor.¹ Oxidation at C-11 of fluprostenol yields 11-keto fluprostenol. 11-keto Fluprostenol exhibits moderate binding to the CRTH2/DP₂ receptor compared to PGD₂ and essentially no activity at the DP₁ receptor.²

References

1. Dukes, M., Russell, W., and Walpole, A.L. Potent luteolytic agents related to prostaglandin F_{2α}. *Nature* **250(464)**, 330-331 (1974).
2. Monneret, G., Cossette, C., Gravel, S., *et al.* 15R-methyl-prostaglandin D₂ is a potent and selective CRTH2/DP₂ receptor agonist in human eosinophils. *J. Pharmacol. Exp. Ther.* **304(1)**, 349-355 (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, 08/15/2022

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