

PRODUCT INFORMATION



Fluprostenol-d₄ Item No. 316767

Formal Name: 9 α ,11 α ,15R-trihydroxy-16-(3-(trifluoromethyl)phenoxy)-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-oic-3,3,4,4-d₄ acid

Synonym: 16-m-trifluoromethylphenoxy tetranor Prostaglandin F_{2 α} -d₄

MF: C₂₃H₂₅D₄F₃O₆

FW: 462.5

Chemical Purity: $\geq 98\%$ (Fluprostenol)

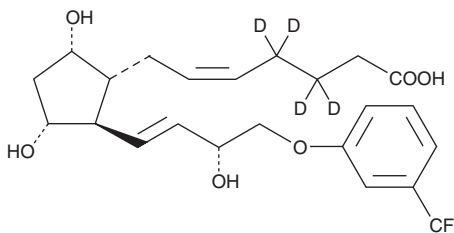
Deuterium Incorporation: $\geq 99\%$ deuterated forms (d₁-d₄); $\leq 1\%$ d₀

UV/Vis.: λ_{max} : 222, 280 nm

Supplied as: A 500 $\mu\text{g/ml}$ solution in methyl acetate

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

Fluprostenol-d₄ is intended for use as an internal standard for the quantification of Fluprostenol (Item No. 16768) 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-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-d₄ in these solvents is approximately 100 mg/ml.

Description

Fluprostenol-d₄ contains four deuterium atoms at the 3, 3', 4, and 4' positions. Fluprostenol is a metabolically stable analog of prostaglandin F_{2 α} (PGF_{2 α}) with potent FP receptor agonist activity.^{1,2} It inhibits PGF_{2 α} binding to human and rat FP receptors with IC₅₀ values of 3.5 and 7.5 nM, respectively.^{1,2} Fluprostenol is a much more potent luteolytic agent than PGF_{2 α} in rats with a minimum fully effective dose of 270 g/kg to terminate pregnancy.³ It is also an effective inhibitor of rat adipose precursor differentiation in primary cultures with an IC₅₀ value of 3-10 $\times 10^{-11}$ M.⁴

References

1. Dukes, M., Russell, W., and Walpole, A.L. Potent luteolytic agents related to prostaglandin F_{2 α} . *Nature* **250** (464), 330-331 (1974).
2. Lake, S., Gullberg, H., Wahlqvist, J., et al. Cloning of the rat and human prostaglandin F_{2 α} receptors and the expression of the rat prostaglandin F_{2 α} receptor. *FEBS Lett.* **355**(3), 317-325 (1994).
3. Abramovitz, M., Boie, Y., Nguyen, T., et al. Cloning and expression of a cDNA for the human prostanoid FP receptor. *J. Biol. Chem.* **269**(4), 2632-2636 (1994).
4. Serrero, G. and Lepak, N.M. Prostaglandin F_{2 α} receptor (FP receptor) agonists are potent adipose differentiation inhibitors for primary culture of adipocyte precursors in defined medium. *Biochem. Biophys. Res. Commun.* **233**(1), 200-202 (1997).

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, 07/13/2026

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