

PRODUCT INFORMATION



16-(3-fluorophenoxy) tetranor Prostaglandin F_{2α} methyl amide

Item No. 26372

Formal Name: (Z)-7-((1R,2R,3R,5S)-2-((R,E)-4-(3-fluorophenoxy)-3-hydroxybut-1-en-1-yl)-3,5-dihydroxycyclopentyl)-N-methylhept-5-enamide

Synonym: 16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide

MF: C₂₃H₃₂FNO₅

FW: 421.5

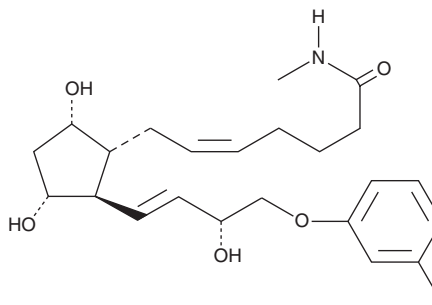
Purity: ≥98%

UV/Vis.: λ_{max}: 220 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

16-(3-fluorophenoxy) tetranor Prostaglandin F_{2α} methyl amide (16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide) 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 16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide in these solvents is approximately 100 mg/ml.

16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide should be diluted with the aqueous buffer of choice. The solubility of 16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

16-(3-fluorophenoxy) tetranor PGF_{2α} methyl amide is a derivative of (±)-cloprostenol (Item No. 16764), a PGF_{2α} analog.

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, 01/31/2024

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