

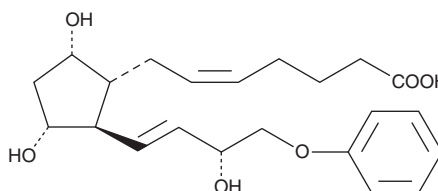
PRODUCT INFORMATION



16-phenoxy tetranor Prostaglandin F_{2α}

Item No. 16760

CAS Registry No.: 51705-19-2
Formal Name: (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(1E,3R)-3-hydroxy-4-phenoxy-1-buten-1-yl]cyclopentyl]-5-heptenoic acid
Synonym: 16-phenoxy tetranor PGF_{2α}
MF: C₂₂H₃₀O₆
FW: 390.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219 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

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

16-phenoxy tetranor PGF_{2α} is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 16-phenoxy tetranor PGF_{2α} should be diluted with the aqueous buffer of choice. The solubility of 16-phenoxy tetranor PGF_{2α} in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

16-phenoxy tetranor PGF_{2α} is a metabolically stable analog of PGF_{2α}. It binds to the FP receptor on ovine luteal cells with much greater affinity (440%) than PGF_{2α}.¹

Reference

1. Balapure, A.K., Rexroad, C.E., Jr., Kawada, K., *et al.* Structural requirements for prostaglandin analog interaction with the ovine corpus luteum prostaglandin F_{2α} receptor. *Biochem. Pharmacol.* **38(14)**, 2375-2381 (1989).

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, 03/15/2024

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