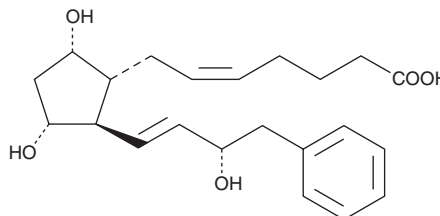


PRODUCT INFORMATION



16-phenyl tetranor Prostaglandin F_{2α} Item No. 16770

CAS Registry No.: 38315-48-9
Formal Name: (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[[1E,3S)-3-hydroxy-4-phenyl-1-buten-1-yl]cyclopentyl]-5-heptenoic acid
Synonym: 16-phenyl tetranor PGF_{2α}
MF: C₂₂H₃₀O₅
FW: 374.5
Purity: ≥99%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

16-phenyl tetranor Prostaglandin F_{2α} (16-phenyl tetranor PGF_{2α}) is supplied as a crystalline solid. A stock solution may be made by dissolving the 16-phenyl tetranor PGF_{2α} in the solvent of choice, which should be purged with an inert gas. 16-phenyl tetranor PGF_{2α} is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 16-phenyl tetranor PGF_{2α} in these solvents is approximately 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 16-phenyl tetranor PGF_{2α} can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 16-phenyl 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-phenyl tetranor PGF_{2α} is a metabolically stable analog of PGF_{2α}. The affinity of 16-phenyl tetranor PGF_{2α} at the FP receptor of ovine luteal cells is poor (8.7%) compared to 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, 02/09/2024

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