

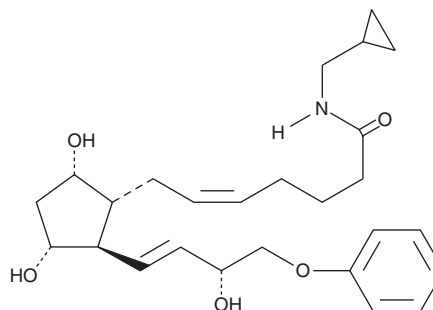
PRODUCT INFORMATION



16-phenoxy tetranor Prostaglandin F_{2α} cyclopropyl methyl amide

Item No. 10010809

CAS Registry No.: 1138395-09-1
Formal Name: 95Z)-N-(cyclopropylmethyl)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(1E,3R)-3-hydroxy-4-phenoxy-1-buten-1-yl]cyclopentyl]-5-heptenamide
Synonym: 16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide
MF: C₂₆H₃₇NO₅
FW: 443.6
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 220, 271 nm
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-phenoxy tetranor Prostaglandin F_{2α} cyclopropyl methyl amide (16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide) is supplied as a crystalline solid. A stock solution may be made by dissolving the 16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide in the solvent of choice, which should be purged with an inert gas. 16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide 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-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide in PBS (pH 7.2) is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

16-phenoxy tetranor PGF_{2α} cyclopropyl methyl amide is an amidated form of 16-phenoxy tetranor PGF_{2α} (Item No. 16760).

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, 07/21/2026

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