

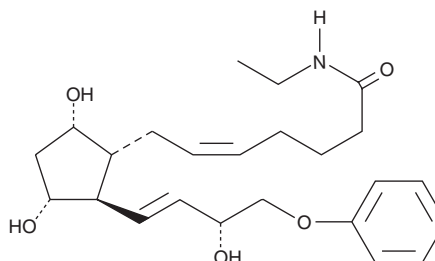
PRODUCT INFORMATION



16-phenoxy Prostaglandin F_{2α} ethyl amide

Item No. 10009875

CAS Registry No.: 951319-59-8
Formal Name: N-ethyl-9α,11α,15S-trihydroxy-16-phenoxy-prosta-5Z,13E-dien-1-amide
Synonym: 16-phenoxy PGF_{2α} ethyl amide
MF: C₂₄H₃₅NO₅
FW: 417.5
Purity: ≥98%
UV/Vis.: λ_{max}: 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 Prostaglandin F_{2α} ethyl amide (16-phenoxy PGF_{2α} ethyl amide) is supplied as a crystalline solid. A stock solution may be made by dissolving the 16-phenoxy PGF_{2α} ethyl amide in the solvent of choice, which should be purged with an inert gas. 16-phenoxy PGF_{2α} ethyl amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 16-phenoxy PGF_{2α} ethyl amide in ethanol is approximately 25 mg/ml and approximately 50 mg/ml in DMSO and DMF.

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 PGF_{2α} ethyl amide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 16-phenoxy PGF_{2α} ethyl amide in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGF_{2α} acting through the FP receptor, causes smooth muscle contraction and exhibits potent luteolytic activity.¹⁻³ 16-phenoxy 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}.⁴ Ethyl amides of PGs serve as prodrugs, as they are hydrolyzed in certain tissues to generate the bioactive free acid.

References

1. Samuelsson, B., Goldyne, M., Granström, E., *et al.* Prostaglandins and thromboxanes. *Annu. Rev. Biochem.* **47**, 997-1029 (1978).
2. Speroff, L. and Ramwell, P.W. Prostaglandins in reproductive physiology. *Am. J. Obstet. Gynecol.* **107**, 1111-1130 (1970).
3. Crankshaw, D.J. and Gaspar, V. Pharmacological characterization *in vitro* of prostanoid receptors in the myometrium of nonpregnant ewes. *J. Reprod. Fertil.* **103**, 55-61 (1995).
4. 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**, 2375-2381 (1989).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY USE - 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, 01/25/2024

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