

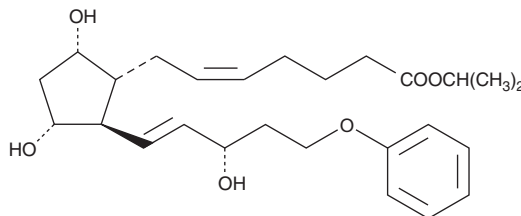
PRODUCT INFORMATION



17-phenoxy trinor Prostaglandin F_{2α} isopropyl ester

Item No. 10010723

Formal Name: 9α,11α,15S-trihydroxy-17-phenoxy-prosta-5Z,13E-dien-1-oic acid, isopropyl ester
Synonym: 17-phenoxy-PGF_{2α} isopropyl ester
MF: C₂₆H₃₈O₆
FW: 446.6
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 221, 272 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

17-phenoxy trinor Prostaglandin F_{2α} isopropyl ester 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 17-phenoxy trinor prostaglandin F_{2α} isopropyl ester in these solvents is approximately 100 mg/ml.

17-phenoxy trinor Prostaglandin F_{2α} isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 17-phenoxy trinor prostaglandin F_{2α} isopropyl ester should be diluted with the aqueous buffer of choice. The solubility of 17-phenoxy trinor prostaglandin F_{2α} isopropyl ester in PBS (pH 7.2) is approximately 100 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

The actions of many clinical F-series prostaglandins (PGs), including those used for estrous synchronization and reduction of intraocular pressure (IOP), are mediated through the PGF_{2α} (FP) receptor.¹⁻³ 17-phenoxy PGF_{2α} isopropyl ester is an analog of PGF_{2α} containing a 17-phenoxy group on the lower side chain and an isopropyl ester at the C-1 position. There are no published reports on the biological activity of 17-phenoxy PGF_{2α} isopropyl ester.

References

1. Narumiya, S., Sugimoto, Y., and Ushikubi, F. Prostanoid receptors: Structures, properties, and functions. *Physiol. Rev.* **79**(4), 1193-1226 (1999).
2. Resul, B., Stjernschantz, J., Selén, G., et al. Structure-activity relationships and receptor profiles of some ocular hypotensive prostanoids. *Surv. Ophthalmol.* **41**(Suppl. 2), S47-S52 (1997).
3. Clark, A.F. and Yorio, T. Ophthalmic drug discovery. *Nature* **2**(6), 448-459 (2003).

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, 04/12/2024

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