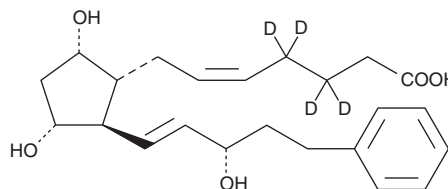


PRODUCT INFORMATION



17-phenyl trinor Prostaglandin F_{2α}-d₄ Item No. 316810

CAS Registry No.: 58976-50-4
Formal Name: 9α,11α,15S-trihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-oic-3,3,4,4-d₄ acid
Synonyms: Bimatoprost (free acid)-d₄, 17-phenyl trinor PGF_{2α}-d₄
MF: C₂₃H₂₈D₄O₅
FW: 392.5
Chemical Purity: ≥98% (17-phenyl trinor PGF_{2α})
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
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

17-phenyl trinor Prostaglandin F_{2α}-d₄ (17-phenyl trinor PGF_{2α}-d₄) is intended for use as an internal standard for the quantification of 17-phenyl trinor PGF_{2α} by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

17-phenyl trinor PGF_{2α}-d₄ 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 (DMF) purged with an inert gas can be used. The solubility of 17-phenyl trinor PGF_{2α}-d₄ in ethanol is approximately 30 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

17-phenyl trinor PGF_{2α} is a metabolically stable analog of PGF_{2α} and is a potent agonist for the FP receptor. It binds to the FP receptor on ovine luteal cells with a relative potency of 756% compared to that of PGF_{2α}.¹ At the rat recombinant FP receptor expressed in CHO cells 17-phenyl trinor PGF_{2α} inhibits PGF_{2α} binding with a K_i of 1.1 nM.² The isopropyl ester of 17-phenyl trinor PGF_{2α}-d₄ is slightly better than PGF_{2α} isopropyl ester in reducing the intraocular pressure in the cat eye without any irritation.³

References

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).
2. Lake, S., Gullberg, H., Wahlqvist, J., *et al.* Cloning of the rat and human prostaglandin F_{2α} receptors and the expression of the rat prostaglandin F_{2α} receptor. *FEBS Lett.* **355(3)**, 317-325 (1994).
3. Stjernschantz, J. and Resul, B. Phenyl substituted prostaglandin analogs for glaucoma treatment. *Drugs Future* **17(8)**, 691-704 (1992).

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, 11/09/2023

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