

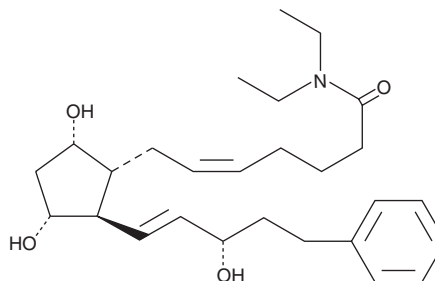
PRODUCT INFORMATION



17-phenyl trinor Prostaglandin F_{2α} diethyl amide

Item No. 16823

CAS Registry No.: 1176637-26-5
Formal Name: N-diethyl-9α,11α,15S-trihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-amide
Synonyms: Bimatoprost diethyl amide, 17-phenyl trinor PGF_{2α} diethyl amide, 17-p-PGF_{2α}-NEt₂
MF: C₂₇H₄₁NO₄
FW: 443.6
Purity: ≥98%
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} diethyl amide (17-phenyl trinor PGF_{2α} diethyl amide) 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} diethyl amide in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Description

17-phenyl trinor PGF_{2α} diethyl amide is an analog of PGF_{2α} in which the C-1 carboxyl group has been modified to an N-diethyl amide. PG esters have been shown to have ocular hypotensive activity.¹ PG N-ethyl amides were recently introduced as alternative PG hypotensive prodrugs.² Studies have shown that bovine and human corneal tissue converts the N-ethyl amides of various PGs to the free acids with a conversion efficiency of about 2.5 μg/g corneal tissue/hr.³ However, dialkyl amides such as 17-phenyl trinor PGF_{2α} diethyl amide are inert to corneal amidase activity, and are not converted in any detectable amount to the corresponding free acids. These compounds may therefore be useful tools in elucidating the claim that PG amides have intrinsic intraocular hypotensive activity.

References

1. Bito, L.Z. Comparison of the ocular hypotensive efficacy of eicosanoids and related compounds. *Exp. Eye Res.* **38(2)**, 181-184 (1984).
2. Woodward, D.F., Krauss, A.H., Chen, J., et al. The pharmacology of bimatoprost (Lumigan™). *Surv. Ophthalmol.* **45(Suppl. 4)**, S337-S345 (2001).
3. Maxey, K.M., Johnson, J., and LaBrecque, J. The hydrolysis of bimatoprost in corneal tissue generates a potent prostanoid FP receptor agonist. *Surv. Ophthalmol.* **47(Suppl. 1)**, S34-S40 (2002).

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, 03/15/2024

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