

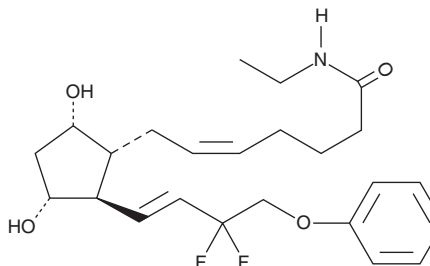
PRODUCT INFORMATION



Tafuprost ethyl amide

Item No. 9000843

CAS Registry No.: 1185851-52-8
Formal Name: N-ethyl-9 α ,11 α -dihydroxy-15,15-difluoro-16-phenoxy-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-amide
Synonyms: Taf EA, Tafuprost EA
MF: C₂₄H₃₃F₂NO₄
FW: 437.5
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Tafuprost ethyl amide is supplied as a solution in ethanol. To change the solvent, simply evaporate the tafuprost ethyl amide under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of tafuprost ethyl amide in these solvents is approximately 30 mg/ml.

Description

Tafuprost ethyl amide is a derivative of the tafuprost (free acid) prodrug tafuprost (Item No. 10005440).¹ Formulations containing tafuprost ethyl amide have been used in the treatment of eyelash hypotrichosis and in cosmetics.

Reference

1. Krupa, M., Chodyński, M., Ostaszewski, A., *et al.* A novel convergent synthesis of the potent antiglaucoma agent tafuprost. *Molecules* **22**(2), 217 (2017).

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. It is the responsibility of the purchaser to determine suitability for other applications.

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, 08/22/2023

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