

PRODUCT INFORMATION



Latanoprost ethyl amide-d₄ Item No. 316822

Formal Name: N-ethyl-9 α ,11 α ,15S-trihydroxy-17-phenyl-18,19,20-trinor-prost-5Z-en-1-
amide-3,3,4,4-d₄

Synonyms: Lat-NEt-d₄

MF: C₂₅H₃₄D₄NO₄

FW: 420.6

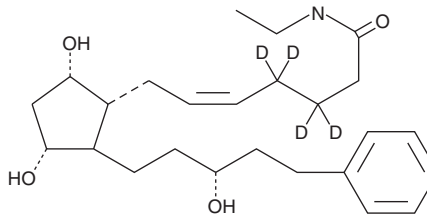
Chemical Purity: $\geq 98\%$ (Latanoprost ethyl amide)

Deuterium Incorporation: $\geq 99\%$ deuterated forms (d₁-d₄); $\leq 1\%$ d₀

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

Latanoprost ethyl amide-d₄ (Lat-NEt-d₄) is intended for use as an internal standard for the quantification of Lat-NEt 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).

Lat-NEt-d₄ 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 Lat-NEt-d₄ in these solvents is approximately 50 and 30 mg/ml, respectively.

Lat-NEt-d₄ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of Lat-NEt-d₄ should be diluted with the aqueous buffer of choice. The solubility of Lat-NEt-d₄ in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lat-NEt-d₄ is intended for use as an internal standard for the quantification of Lat-NEt by GC- or LC-mass spectrometry. Lat-NEt is an F-series prostaglandin (PG) analog in which the C-1 carboxyl group has been modified to an N-ethyl amide. PG esters have been shown to have ocular hypotensive activity.¹ PG N-ethyl amides were recently introduced as alternative PG ocular hypotensive prodrugs.² Although it has been claimed that PG ethyl amides are not converted to the free acids *in vivo*,² studies in our laboratories have shown that bovine and human corneal tissue converts the N-ethyl amides of various PGs to the free acids with a conversion rate of about 2.5 $\mu\text{g/g}$ corneal tissue/hr.³ Lat-NEt would be expected to show the typical intraocular effects of Lat free acid, but with the much slower hydrolysis pharmacokinetics of the PG N-amides.

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.-P., Chen, J., *et al.* The pharmacology of Bimatoprost (LumiganTM). *Surv. Ophthalmol.* **45 (Suppl. 4)**, S337-S345 (2001).
3. Maxey, K.M., Johnson, J., Camras, C.B., *et al.* The hydrolysis of bimatoprost in corneal tissue generates a potent prostanoid FP receptor agonist. *Surv. Ophthalmol.* **47(4)**, 34-40 (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, 04/02/2024

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