

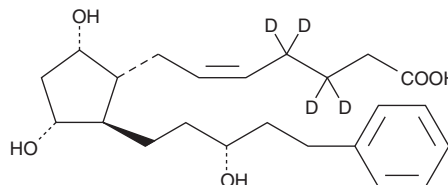
# PRODUCT INFORMATION



## Latanoprost (free acid)-d<sub>4</sub>

Item No. 316811

**CAS Registry No.:** 1224443-47-3  
**Formal Name:** 9 $\alpha$ ,11 $\alpha$ ,15R-trihydroxy-17-phenyl-18,19,20-trinor-prost-5Z-en-1-oic-3,3,4,4-d<sub>4</sub> acid  
**Synonyms:** 17-phenyl-13,14-dihydro trinor Prostaglandin F<sub>2 $\alpha$</sub> -d<sub>4</sub>, PhXA 85-d<sub>4</sub>, Lat-FA-d<sub>4</sub>  
**MF:** C<sub>23</sub>H<sub>30</sub>D<sub>4</sub>O<sub>5</sub>  
**FW:** 394.5  
**Chemical Purity:**  $\geq$ 98% (Latanoprost (free acid))  
**Deuterium Incorporation:**  $\geq$ 99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>);  $\leq$ 1% d<sub>0</sub>  
**Supplied as:** A solution in methyl acetate  
**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

Latanoprost (free acid)-d<sub>4</sub> is intended for use as an internal standard for the quantification of latanoprost (free acid) (Item No. 16811) 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).

Latanoprost (free acid)-d<sub>4</sub> 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 latanoprost (free acid)-d<sub>4</sub> in ethanol is approximately 100 mg/ml and approximately 50 mg/ml in DMSO and DMF.

### Description

Latanoprost (free acid)-d<sub>4</sub> is intended for use as an internal standard for the quantification of latanoprost (free acid) (Item No. 16811) by GC- or LC-MS. Latanoprost (free acid) is a derivative of prostaglandin F<sub>2 $\alpha$</sub>  (PGF<sub>2 $\alpha$</sub> ; Item Nos. 16010 | 16020), an FP receptor agonist, and an active metabolite of the prodrug latanoprost (Item No. 16812).<sup>1</sup> It selectively binds to the FP receptor (K<sub>i</sub> = 0.098  $\mu$ M) over the EP<sub>1</sub>, EP<sub>2</sub>, EP<sub>3</sub>, and EP<sub>4</sub> receptors (K<sub>s</sub> = 2.06, 39.667, 7.519, and 75  $\mu$ M, respectively), as well as the DP, IP, and TP receptors (K<sub>s</sub> =  $\geq$ 20,  $\geq$ 90, and  $\geq$ 60  $\mu$ M, respectively). It induces phosphoinositide turnover in isolated human ciliary muscle and human trabecular meshwork cells, mouse NIH3T3 fibroblasts, and rat A7r5 vascular smooth muscle cells, which all endogenously express FP receptors (EC<sub>50</sub>s = 124, 35, 32, and 35 nM, respectively), as well as HEK293 cells expressing human ocular FP receptors (EC<sub>50</sub> = 45.7 nM). It also increases intracellular calcium levels in hEP<sub>1</sub>-5/293-AEQ17 cells expressing EP<sub>1</sub> receptors in an aequorin-based calcium assay (EC<sub>50</sub> = 119 nM).<sup>2</sup> Latanoprost (free acid) induces contraction of isolated cat iris sphincter smooth muscle (EC<sub>50</sub> = 29.9 nM).<sup>3</sup>

#### 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

# PRODUCT INFORMATION



## References

---

1. Sharif, N.A., Kelly, C.R., Crider, J.Y., *et al.* Ocular hypotensive FP prostaglandin (PG) analogs: PG receptor subtype binding affinities and selectivities, and agonist potencies at FP and other PG receptors in cultured cells. *J. Ocul. Pharmacol. Ther.* **19(6)**, 501-515 (2003).
2. Gurney, M.E., Nugent, R.A., Mo, X., *et al.* Design and synthesis of selective phosphodiesterase 4D (PDE4D) allosteric inhibitors for the treatment of Fragile X Syndrome and other brain disorders. *J. Med. Chem.* **62(10)**, 4884-4901 (2019).
3. Sharif, N.A., Kaddour-Djebbar, I., and Abdel-Latif, A.A. Cat iris sphincter smooth-muscle contraction: Comparison of FP-class prostaglandin analog agonist activities. *J. Ocul. Pharmacol. Ther.* **24(2)**, 152-163 (2008).

## 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