

# PRODUCT INFORMATION

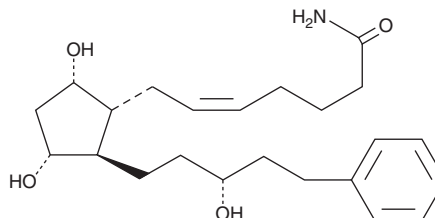


## Latanoprost amide

Item No. 25245

**CAS Registry No.:** 919281-21-3  
**Formal Name:** (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(3R)-3-hydroxy-5-phenylpentyl]cyclopentyl]-5-heptenamide

**MF:** C<sub>23</sub>H<sub>35</sub>NO<sub>4</sub>  
**FW:** 389.5  
**Purity:** ≥98%  
**Supplied as:** An oil  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

Latanoprost amide is supplied as an oil. A stock solution may be made by dissolving the latanoprost amide in the solvent of choice, which should be purged with an inert gas. Latanoprost amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of latanoprost amide in these solvents is approximately 100 mg/ml in ethanol and DMF and approximately 50 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of latanoprost amide can be prepared by directly dissolving the oil in aqueous buffers. The solubility of latanoprost amide in PBS, pH 7.2, is approximately 0.05 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Latanoprost amide is a derivative of latanoprost (Item No. 16812).

#### 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, 03/24/2021

#### 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