

PRODUCT INFORMATION



Latanoprost dimethyl amide

Item No. 25247

Formal Name: (Z)-7-((1R,2R,3R,5S)-3,5-dihydroxy-2-((R)-3-hydroxy-5-phenylpentyl)cyclopentyl)-N,N-dimethylhept-5-enamide

MF: C₂₅H₃₉NO₄

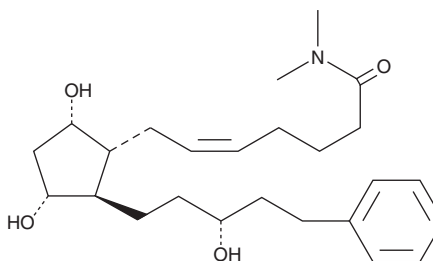
FW: 417.6

Purity: ≥98%

Supplied as: A neat 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 dimethyl amide is supplied as a neat oil. A stock solution may be made by dissolving the latanoprost dimethyl amide in the solvent of choice, which should be purged with an inert gas. Latanoprost dimethyl amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of latanoprost dimethyl amide in ethanol and DMF is approximately 100 mg/ml 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 dimethyl amide can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of latanoprost dimethyl 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 dimethyl amide is a derivative of latanoprost (Item No. 16812) and an isomer of latanoprost ethyl amide (Item No. 16822).

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, 05/21/2020

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