

# PRODUCT INFORMATION



## 9-keto Latanoprost

Item No. 9000350

**CAS Registry No.:** 141074-06-8  
**Formal Name:** (5Z)-7-[(1R,2R,3R)-3-hydroxy-2-[(3R)-3-hydroxy-5-phenylpentyl]-5-oxocyclopentyl]-5-heptenoic acid, 1-methylethyl ester

**MF:** C<sub>26</sub>H<sub>38</sub>O<sub>5</sub>

**FW:** 430.6

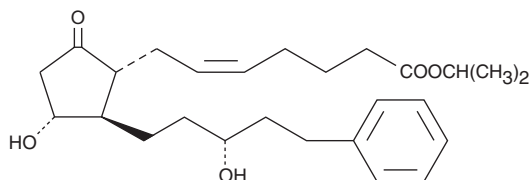
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 213, 261, 269 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

9-keto Latanoprost is supplied as a crystalline solid. A stock solution may be made by dissolving the 9-keto latanoprost in the solvent of choice, which should be purged with an inert gas. 9-keto Latanoprost is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 9-keto latanoprost in these solvents is approximately 100 mg/ml.

9-keto Latanoprost is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 9-keto latanoprost should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 9-keto Latanoprost has a solubility of approximately 80 µg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

9-keto Latanoprost is a derivative of the FP receptor agonist 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, 04/10/2024

#### 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