

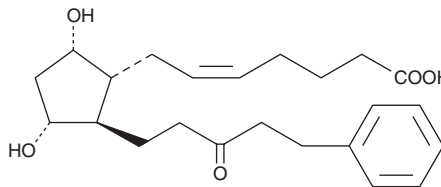
# PRODUCT INFORMATION



## 15-keto Latanoprost (free acid)

Item No. 16815

**CAS Registry No.:** 369585-22-8  
**Formal Name:** (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-(3-oxo-5-phenylpentyl)cyclopentyl]-5-heptenoic acid  
**MF:** C<sub>23</sub>H<sub>32</sub>O<sub>5</sub>  
**FW:** 388.5  
**Purity:** ≥95%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

15-keto Latanoprost (free acid) 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 15-keto latanoprost (free acid) in ethanol and DMF is approximately 50 mg/ml and approximately 33 mg/ml in DMSO.

15-keto Latanoprost (free acid) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15-keto latanoprost (free acid) should be diluted with the aqueous buffer of choice. 15-keto Latanoprost (free acid) has a solubility of 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

15-keto Latanoprost (free acid) is a metabolite of latanoprost (Item No. 16812).<sup>1</sup> It is formed from latanoprost (free acid) via dehydrogenation at position 15 by 15-hydroxyprostaglandin dehydrogenase in monkey cornea.

### Reference

1. Fujimori, K., Okada, T., and Urade, Y. Expression of NADP<sup>+</sup>-dependent 15-hydroxyprostaglandin dehydrogenase mRNA in monkey ocular tissues and characterization of its recombinant enzyme. *J. Biochem.* **131**(3), 383-389 (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, 03/15/2024

#### 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