

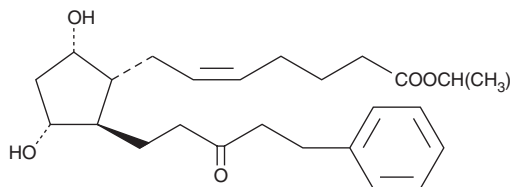
PRODUCT INFORMATION



15-keto Latanoprost

Item No. 16816

CAS Registry No.: 135646-98-9
Formal Name: 9 α ,11 α -dihydroxy-15-oxo-17-phenyl-18,19,20-trinor-prost-5Z-en-1-oic acid, isopropyl ester
MF: C₂₆H₃₈O₅
FW: 430.6
Purity: \geq 98%
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

15-keto Latanoprost 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 purged with an inert gas can be used. The solubility of 15-keto latanoprost in these solvents is approximately 100 mg/ml.

15-keto Latanoprost is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15-keto latanoprost should be diluted with the aqueous buffer of choice. 15-keto Latanoprost has a solubility of 0.08 mg/ml in a 1:12 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 is an active metabolite of the FP receptor agonist latanoprost (Item No. 16812).^{1,2} Intraocular administration of 15-keto latanoprost (1 μ g/eye) reduces intraocular pressure in cynomolgus monkeys.² It also reduces pupillary diameter in cats when administered topically at a dose of 5 μ g/eye. 15-keto Latanoprost is a potential pharmaceutical impurity found in commercial preparations of latanoprost.

References

1. Fujimori, K., Okada, T., and Urade, Y. Expression of NADP⁺-dependent 15-hydroxyprostaglandin dehydrogenase mRNA in monkey ocular tissues and characterization of its recombinant enzyme. *J. Biochem.* **131**(3), 383-389 (2002).
2. Stjernschantz, J. and Resul, B. Phenyl substituted prostaglandin analogs for glaucoma treatment. *Drugs Future* **17**(8), 691-704 (1992).

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