

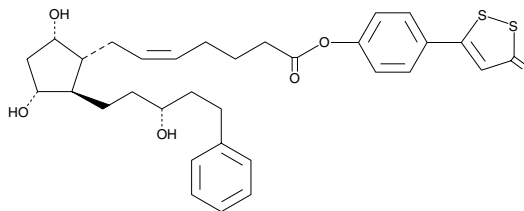
Product Information



ACS 67

Item No. 13619

CAS Registry No.: 1088434-86-9
Formal Name: 7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(3R)-3-hydroxy-5-phenylpentyl]cyclopentyl]-4-(3-thioxo-3H-1,2-dithiol-5-yl)phenyl ester, 5Z-heptenoic acid
MF: C₃₂H₃₈O₅S₃
FW: 598.8
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that ACS 67 be stored as supplied at -20°C. It should be stable for at least two years.

ACS 67 is supplied as a crystalline solid. A stock solution may be made by dissolving the ACS 67 in the solvent of choice. ACS 67 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ACS 67 in these solvents is approximately 2, 20, and 30 mg/ml, respectively.

If aqueous stock solutions are required for biological experiments, they can best be prepared by dissolving the ACS 67 with aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Latanoprost is an F-series prostaglandin (PG) analog which has been approved for use as an ocular hypotensive drug for the treatment of glaucoma.¹ Latanoprost is an isopropyl ester, a prodrug form which is converted to latanoprost (free acid) by endogenous esterase enzymes. The free acid form is 200 times more potent than latanoprost as an FP receptor ligand for the human recombinant FP receptor.² ACS 67 is an analog of latanoprost (free acid) that contains a hydrogen sulfide releasing component conjugated to the latanoprost carboxyl group.³ In glaucomatous pigmented rabbits, ACS 67 reduced intraocular pressure (IOP) more rapidly and to a greater extent than latanoprost (15 *versus* 25.5 mm Hg at four hours) at a dose of 0.005% for each compound.³ ACS 67 also increased the levels of reduced glutathione and cGMP in the aqueous humor of glaucomatous pigmented rabbits better than latanoprost.³

References

1. Stjernschantz, J. and Resul, B. Phenyl substituted prostaglandin analogs for glaucoma treatment. *Drugs of the Future* **17**, 691-704 (1992).
2. Abramovitz, M., Adam, M., Boie, Y., *et al.* The utilization of recombinant prostanoid receptors to determine the affinities and selectivities of prostaglandins and related analogs. *Biochim. Biophys. Acta* **1483**, 285-293 (2000).
3. Perrino, E., Uliva, C., Lanzi, C., *et al.* New prostaglandin derivative for glaucoma treatment. *Bioorg. Medicinal Chem. Letters* **19**, 1639-1642 (2009).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13619

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/02/2011