

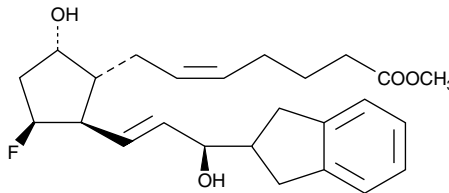
Product Information



AL 8810 methyl ester

Item No. 10008370

CAS Registry No.: 1176541-11-9
Formal Name: 9 α ,15R-dihydroxy-11 β -fluoro-15-(2,3-dihydro-1H-inden-2-yl)-16,17,18,19,20-pentanoic acid, methyl ester
MF: C₂₅H₃₃FO₄
FW: 416.5
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that AL 8810 methyl ester be stored as supplied at -20°C. It should be stable for at least two years.

AL 8810 methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the AL 8810 methyl ester in an organic solvent purged with an inert gas. AL 8810 methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AL 8810 methyl ester in these solvents is approximately 50, 25, and 30 mg/ml, respectively.

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 AL 8810 methyl ester can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of AL 8810 methyl ester in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

AL 8810 is an 11 β -fluoro analog of prostaglandin F_{2 α} (PGF_{2 α}) which acts as a potent and selective antagonist at the FP receptor.¹ AL 8810 methyl ester is a lipid soluble, esterified prodrug form of AL 8810 analogous to the commonly used therapeutic intraocular PG compounds such as latanoprost² and travoprost.³ The pharmacology of AL 8810 methyl ester has not been published.

References

1. Griffen, B.W., Klimko, P., Crider, J.Y., *et al.* AL-8810: A novel prostaglandin F_{2 α} analog with selective antagonist effects at the prostaglandin F_{2 α} (FP) receptor. *J. Phar. Exp. Ther.* **290**, 1278-1284 (1999).
2. Camras, C.B., Alm, A., Watson, P., *et al.* Latanoprost, a prostaglandin analog, for glaucoma therapy. Efficacy and safety after 1 year of treatment in 198 patients. *Ophthalmology* **103**, 1916-1924 (1996).
3. Sorbera, L.A. and Castañer, J. Travoprost. *Drugs of the Future* **25**, 41-45 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008370

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.

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 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, 11/22/2013