

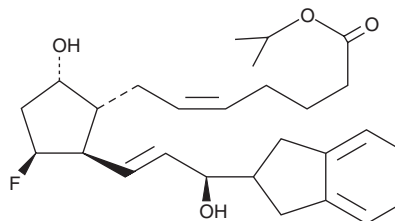
# PRODUCT INFORMATION



## AL 8810 isopropyl ester

Item No. 10113

**CAS Registry No.:** 208114-93-6  
**Formal Name:** 9 $\alpha$ ,15R-dihydroxy-11 $\beta$ -fluoro-15-(2,3-dihydro-1H-inden-2-yl)-16,17,18,19,20-pentanoic acid, isopropyl ester  
**MF:** C<sub>27</sub>H<sub>37</sub>FO<sub>4</sub>  
**FW:** 444.6  
**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 isopropyl ester be stored as supplied at -20°C. It should be stable for at least two years.

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

AL 8810 isopropyl is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AL 8810 isopropyl should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. AL 8810 isopropyl has a solubility of 100  $\mu$ g/ml in approximately 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

AL 8810 (Item No. 16735) is an 11 $\beta$ -fluoro analog of prostaglandin F<sub>2 $\alpha$</sub>  (Item No. 16010) which acts as a potent and selective antagonist at the FP receptor.<sup>1</sup> AL 8810 isopropyl ester is a lipid soluble, esterified prodrug form of AL 8810 analogous to the commonly used therapeutic intraocular prostaglandin compounds such as Latanoprost (Item No. 16812) and Travoprost.<sup>2,3</sup> The pharmacology of AL 8810 isopropyl ester has not been published.

### References

1. Griffen, B.W., Klimko, P., Crider, J.Y., *et al.* AL-8810: A novel prostaglandin F<sub>2 $\alpha$</sub>  analog with selective antagonist effects at the prostaglandin F<sub>2 $\alpha$</sub>  (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).

#### 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, 08/04/2015

#### 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