

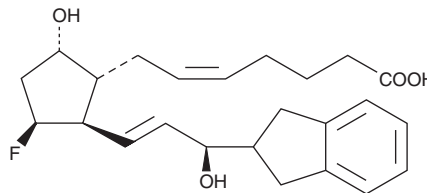
# PRODUCT INFORMATION



## AL 8810

Item No. 16735

**CAS Registry No.:** 246246-19-5  
**Formal Name:** 9 $\alpha$ ,15R-dihydroxy-11 $\beta$ -fluoro-15-(2,3-dihydro-1H-inden-2-yl)-16,17,18,19,20-pentanoic acid  
**MF:** C<sub>24</sub>H<sub>31</sub>O<sub>4</sub>F  
**FW:** 402.5  
**Purity:**  $\geq$ 98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

AL 8810 is supplied as a crystalline solid. A stock solution may be made by dissolving the AL 8810 in the solvent of choice. AL 8810 is miscible in the organic solvent ethanol, which should be purged with an inert gas. AL 8810 is soluble in DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of AL 8810 in these solvents is approximately 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 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of AL 8810 in PBS, pH 7.2, is approximately 0.05 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

AL 8810 is an 11 $\beta$ -fluoro analog of PGF<sub>2 $\alpha$</sub>  which acts as a potent and selective antagonist at the FP receptor.<sup>1</sup> AL 8810 has weak intrinsic agonist activity on FP receptor preparations in the 200-300 nM range. Yet it fully antagonized the activity of the potent FP receptor agonist fluprostenol at this concentration, with EC<sub>50</sub> values of approximately 430 nM. AL 8810 fully antagonized the bimatoprost-induced calcium mobilization in Swiss 3T3 fibroblasts at 100  $\mu$ M, indicating that bimatoprost acts as an FP agonist in this preparation.<sup>2</sup> The K<sub>i</sub> for the inhibition of several potent agonists at the cloned human ciliary body FP receptor is in the range of 1-2  $\mu$ M.<sup>3</sup>

### 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. Pharm. Exp. Ther.* **290**, 1278-1284 (1999).
2. Sharif, N.A., Williams, G.W., and Kelly, C.R. Bimatoprost and its free acid are prostaglandin FP receptor agonists. *Eur. J. Pharmacol.* **432**, 211-213 (2001).
3. Sharif, N.A., Kelly, C.R., and Crider, J.Y. Agonist activity of Bimatoprost, Travoprost, Latanoprost, Unoprostone isopropyl ester and other prostaglandin analogs at the cloned human ciliary body FP prostaglandin receptor. *J. Ocul. Pharmacol.* **18**, 313-324 (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, 02/09/2024

#### 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