

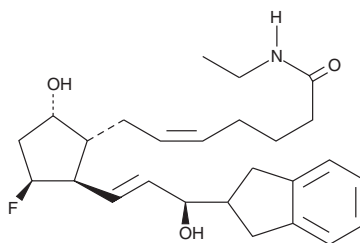
PRODUCT INFORMATION



AL 8810 ethyl amide

Item No. 10010470

Formal Name: 9 α ,15R-dihydroxy-11 β -fluoro-15-92,3-dihydro-1H-inden-2-yl)-16,17,18,19,20-pentanor-prosta-5Z,13E-dion-1-oic acid, ethyl amide
MF: C₂₆H₃₆FNO₃
FW: 429.6
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

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

AL 8810 ethyl amide is supplied as a crystalline solid. A stock solution may be made by dissolving the AL 8810 ethyl amide in an organic solvent purged with an inert gas. AL 8810 ethyl amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AL 8810 ethyl amide in ethanol and DMF is approximately 15 mg/ml and approximately 10 mg/ml in DMSO.

AL 8810 ethyl amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AL 8810 ethyl amide should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. AL 8810 ethyl amide has a solubility of approximately 0.3 mg/ml in a 1:2 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 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 ethyl amide is an analog of AL 8810 in which the C-1 carboxyl group has been modified to an N-ethyl amide. This modification is analogous to the PG N-ethyl amides, as typified by Bimatoprost, that have been introduced as alternative PG ocular hypotensive prodrugs.² In contrast to AL 8810 which contracted the cat iris, AL 8810 ethyl amide showed no contraction activity at concentrations up to 10⁻⁴ M and did not antagonize the activity of PGF_{2 α} -ethanolamide in this system.³

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. Pharmacol. Exp. Ther.* **290**, 1278-1284 (1999).
2. Woodward, D.F., Krauss, A.H.-P., Chen, J., *et al.* The pharmacology of BimatoprostTM (LumiganTM). *Survey of Ophthalmology* **45**, S337-S345 (2001).
3. Woodward, D.F., Krauss, A.H., Wang, J.W., *et al.* Identification of an antagonist that selectively blocks the activity of prostamides (prostaglandin-ethanolamides) in the feline iris. *Br. J. Pharmacol.* **150**, 342-352 (2007).

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, 12/22/2015

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