

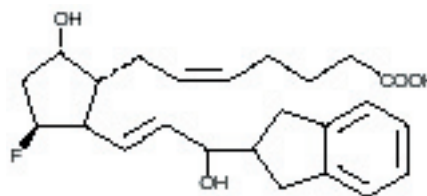
Product Information



(S)-AL 8810

Catalog No. 10023

Formal Name: 9 α ,15S-dihydroxy-11 β -fluoro-15-(2,3-dihydro-1H-inden-2-yl)-16,17,18,19,20-pentanoic-acid
MF: C₂₄H₃₁O₄F
FW: 402.5
Purity: \geq 98%
Stability: \geq 1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that (S)-AL 8810 be stored as supplied at -20°C. It will be stable for at least one year.

(S)-AL 8810 is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-AL 8810 in an organic solvent purged with an inert gas. (S)-AL 8810 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (S)-AL 8810 in these solvents is approximately 25 mg/ml. (S)-AL 8810 will be stable for at least six months in these solvents if stored at -20°C.

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 (S)-AL 8810 can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of (S)-AL 8810 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 (S)-AL 8810 is the C-15 epimer of AL 8810, having the inverse, (S), or 'natural' configuration at C-15 relative to AL 8810, which is 15-(R).

The pharmacology of (S)-AL 8810 has not been published.

Reference

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).

Related Products

Prostaglandin F_{2 α} - Cat. No. 16010 • AL 8810 - Cat. No. 16735

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

Copyright Cayman Chemical Company, 03/16/2006

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