

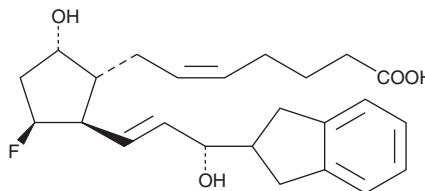
PRODUCT INFORMATION



(S)-AL 8810

Item No. 10023

CAS Registry No.: 1224444-55-6
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 should 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.

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 (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).

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, 07/31/2015

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