

PRODUCT INFORMATION

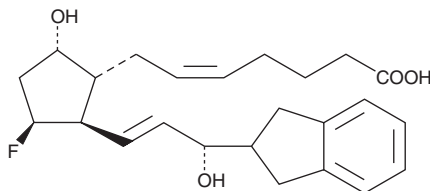


(S)-AL 8810

Item No. 10023

CAS Registry No.: 1224444-55-6
Formal Name: (5Z)-7-[(1R,2R,3S,5S)-2-[(1E,3S)-3-(2,3-dihydro-1H-inden-2-yl)-3-hydroxy-1-propen-1-yl]-3-fluoro-5-hydroxycyclopentyl]-5-heptenoic acid

MF: C₂₄H₃₁O₄F
FW: 402.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(S)-AL 8810 is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-AL 8810 in the solvent of choice, which should be purged with an inert gas. (S)-AL 8810 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (S)-AL 8810 in these solvents is approximately 25 and 30 mg/ml, respectively. (S)-AL 8810 is also miscible in ethanol.

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 solid 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 11β-fluoro analog of PGF_{2α} 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_{2α} analog with selective antagonist effects at the prostaglandin F_{2α} (FP) receptor. *J. Pharmacol. Exp. Ther.* **290(3)**, 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, 04/24/2024

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