

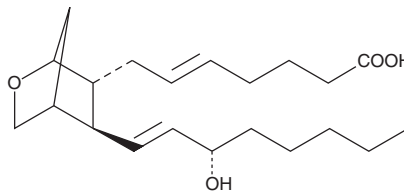
PRODUCT INFORMATION



5-*trans* U-44069

Item No. 16442

CAS Registry No.: 330796-57-1
Formal Name: 9,11-dideoxy-9 α ,11 α -epoxymethano-prosta-5E,13E-dien-1-oic acid
Synonym: 5,6-*trans* U-44069
MF: C₂₁H₃₄O₄
FW: 350.5
Purity: \geq 98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

5-*trans* U-44069 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 5-*trans* U-44069 in these solvents is approximately 100 mg/ml. 5-*trans* U-44069 is 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. If an organic solvent-free solution of 5-*trans* U-44069 is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 5-*trans* U-44069 in PBS (pH 7.2) is at least 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-*trans* U-44069 is the *trans* isomer of the thromboxane receptor agonist U-44069 (Item No. 16440). It inhibits microsomal prostaglandin E₂ synthase (mPGES) when used at a concentration of 10 μ M.¹

Reference

1. Quraishi, O., Mancini, J. A., and Riendeau, D. Inhibition of inducible prostaglandin E₂ synthase by 15-deoxy- $\Delta^{12,14}$ -prostaglandin J₂ and polyunsaturated fatty acids. *Biochem. Pharmacol.* **63**(6), 1183-1189 (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, 06/08/2020

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