

PRODUCT INFORMATION



U-44069 serinol amide

Item No. 40872

Formal Name: (Z)-N-(2,3-dihydroxypropyl)-7-((1S,4R,5S,6R)-5-((S,E)-3-hydroxyoct-1-en-1-yl)-2-oxabicyclo[2.2.1]heptan-6-yl)hept-5-enamide

Synonyms: 9,11-dideoxy-9 α ,11 α -epoxymethano PGF_{2 α} serinol amide, 9,11-dideoxy-9 α ,11 α -epoxymethano Prostaglandin F_{2 α} serinol amide, 9,11-epoxymethano Prostaglandin H₂ serinol amide

MF: C₂₄H₄₁NO₅

FW: 423.6

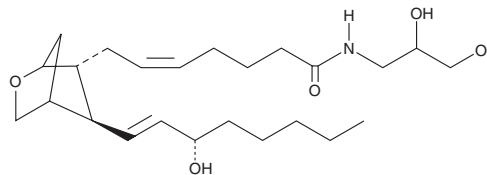
Purity: \geq 95%

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: \geq 1 year

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



Laboratory Procedures

U-44069 serinol amide 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. U-44069 serinol amide is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

U-44069 serinol amide is a derivative of the TP receptor agonist U-44069 (Item No. 16440).

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, 08/06/2024

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