

# PRODUCT INFORMATION



## U-46619 serinol amide

Item No. 40870

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

**Synonyms:** 9,11-dideoxy-9 $\alpha$ ,11 $\alpha$ -methanoepoxy PGF<sub>2 $\alpha$</sub>  serinol amide, 9,11-dideoxy-9 $\alpha$ ,11 $\alpha$ -methanoepoxy Prostaglandin F<sub>2 $\alpha$</sub>  serinol amide

**MF:** C<sub>24</sub>H<sub>41</sub>NO<sub>5</sub>

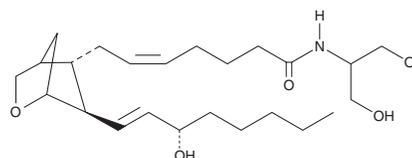
**FW:** 423.6

**Purity:**  $\geq$ 95%

**Supplied as:** A 5 mg/ml solution in methyl acetate

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

U-46619 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. Solvents such as ethanol and DMSO purged with an inert gas can be used. U-46619 serinol amide is soluble ( $\geq$ 10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

### Description

U-46619 serinol amide is a derivative of the TP receptor agonist U-46619 (Item No. 16450).

#### 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, 11/18/2025

#### 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