

PRODUCT INFORMATION



17-phenyl trinor Prostaglandin F_{2α} glycinamide methyl ester

Item No. 25471

Formal Name: methyl ((Z)-7-((1R,2R,3R,5S)-3,5-dihydroxy-2-((S,E)-3-hydroxy-5-phenylpent-1-en-1-yl)cyclopentyl)hept-5-enoyl)glycinate

Synonym: 17-phenyl trinor PGF_{2α} glycinamide methyl ester

MF: C₂₆H₃₇NO₆

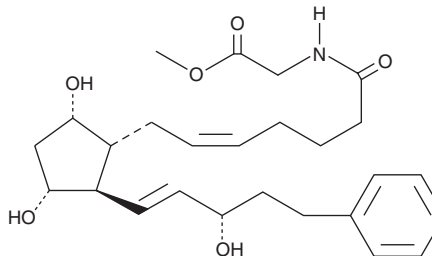
FW: 459.6

Purity: ≥98%

Supplied as: An oil

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

17-phenyl trinor Prostaglandin F_{2α} glycinamide methyl ester is supplied as an oil. A stock solution may be made by dissolving the 17-phenyl trinor prostaglandin F_{2α} glycinamide methyl ester in the solvent of choice, which should be purged with an inert gas. 17-phenyl trinor Prostaglandin F_{2α} glycinamide methyl ester is soluble in organic solvents such as ethanol, acetonitrile, DMSO, and dimethyl formamide (DMF). The solubility of 17-phenyl trinor prostaglandin F_{2α} glycinamide methyl ester in ethanol and acetonitrile is approximately 50 and 3 mg/ml, respectively, and approximately 25 mg/ml in DMSO and DMF.

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 17-phenyl trinor prostaglandin F_{2α} glycinamide methyl ester can be prepared by directly dissolving the oil in aqueous buffers. The solubility of 17-phenyl trinor prostaglandin F_{2α} glycinamide methyl ester in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

17-phenyl trinor Prostaglandin F_{2α} glycinamide methyl ester is a derivative of 17-phenyl trinor prostaglandin F_{2α} ethyl amide (Item No. 16820).

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, 05/21/2020

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