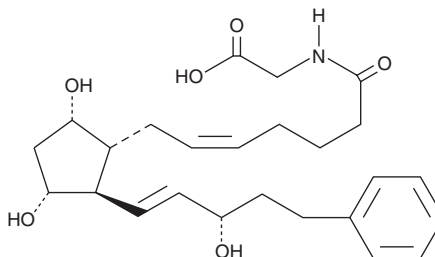


PRODUCT INFORMATION

17-phenyl trinor Prostaglandin F_{2α} glycinamide

Item No. 25470

Formal Name: ((Z)-7-((1R,2R,3R,5S)-3,5-dihydroxy-2-((S,E)-3-hydroxy-5-phenylpent-1-en-1-yl)cyclopentyl)hept-5-enoyl)glycine
Synonym: 17-phenyl trinor PGF_{2α} glycinamide
MF: C₂₅H₃₅NO₆
FW: 445.6
Purity: ≥95%
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 (17-phenyl trinor PGF_{2α} glycinamide) is supplied as an oil. A stock solution may be made by dissolving the 17-phenyl trinor PGF_{2α} glycinamide in the solvent of choice, which should be purged with an inert gas. 17-phenyl trinor PGF_{2α} glycinamide is soluble in organic solvents such as ethanol, acetonitrile, DMSO, and dimethyl formamide (DMF). The solubility of 17-phenyl trinor PGF_{2α} glycinamide 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 PGF_{2α} glycinamide can be prepared by directly dissolving the oil in aqueous buffers. The solubility of 17-phenyl trinor PGF_{2α} glycinamide 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 PGF_{2α} glycinamide is a derivative of 17-phenyl trinor PGF_{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, 05/28/2020

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