

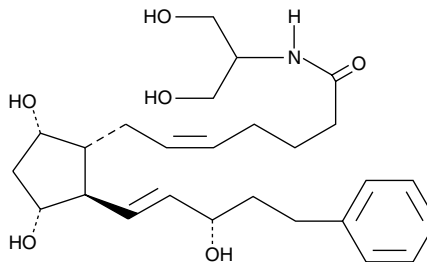
Product Information



17-phenyl trinor Prostaglandin F_{2α} serinol amide

Item No. 10004237

CAS Registry No.: 1175838-84-2
Formal Name: N-[(2-hydroxy-1-hydroxymethyl)ethyl]-9α,11α,15S-trihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-amide
Synonyms: Bimatoprost serinol amide, 17-phenyl trinor PGF_{2α}-SA
MF: C₂₆H₃₉NO₆
FW: 461.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 17-phenyl trinor prostaglandin F_{2α} serinol amide (17-phenyl trinor PGF_{2α} serinol amide) be stored as supplied at -20°C. It should be stable for at least one year.

17-phenyl trinor PGF_{2α} serinol amide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol 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 17-phenyl trinor PGF_{2α} serinol amide in these solvents is at least 10 mg/ml.

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 17-phenyl trinor PGF_{2α} serinol amide is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 17-phenyl trinor PGF_{2α} serinol amide in PBS (pH 7.2) is at least 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

2-Arachidonyl glycerol (2-AG) exhibits cannabinoid (CB) agonist activity at the CB₁ receptor,¹ is an important endogenous monoglyceride species,² and is thus considered to be the natural ligand for the CB₁ receptor. 2-AG can also be metabolized by COX-2 and specific PG synthases to form PG 2-glycerol esters.³ These 2-glycerol esters rapidly equilibrate to form 90:10 mixtures favoring the more stable 1-glycerol ester. 17-phenyl trinor PGF_{2α} serinol amide is a stable analog of PGF_{2α} 2-glycerol ester incorporating an activity-enhancing 17-phenyl substitution. The biological activity of 17-phenyl trinor PGF_{2α} serinol amide has not yet been determined.

References

1. Sugiura, T., Kodaka, T., Kondo, S., *et al. J. Biochem.* **122**, 890-895 (1997).
2. Kondo, S., Kondo, H., Nakane, S., *et al. FEBS Lett.* **429**, 152-156 (1998).
3. Kozak, K.R., Crews, B.C., Morrow, J.D., *et al. J. Biol. Chem.* **277(47)**, 44877-44885 (2002).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10004237

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants **only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/17/2013

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com