

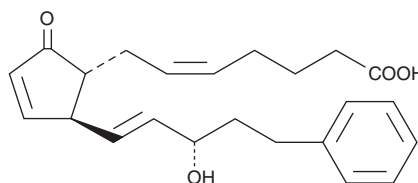
PRODUCT INFORMATION



17-phenyl trinor Prostaglandin A₂

Item No. 10288

CAS Registry No.: 38315-51-4
Formal Name: 9-oxo-15S-hydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,10,13E-trien-1-oic acid
Synonym: 17-phenyl trinor PGA₂
MF: C₂₃H₂₈O₄
FW: 368.5
Purity: ≥97%
UV/Vis.: λ_{max}: 209 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

17-phenyl trinor PGA₂ 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, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 17-phenyl trinor PGA₂ in these solvents is approximately 100 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 PGA₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 17-phenyl trinor PGA₂ in PBS (pH 7.2) is approximately 800 µg/ml. Store aqueous solutions of 17-phenyl trinor PGA₂ on ice and use within 12 hours of preparation. Although the aqueous solutions of 17-phenyl trinor PGA₂ may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Description

17-phenyl trinor PGA₂ is a synthetic prostaglandin analog designed for increased half-life and greater potency. There are no published reports of the biological activity of 17-phenyl trinor PGA₂ in the literature at this time.

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, 03/07/2024

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