

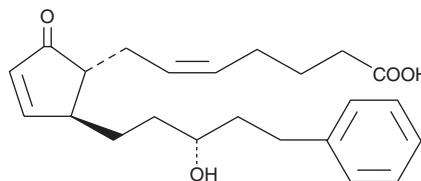
PRODUCT INFORMATION



17-phenyl trinor-13,14-dihydro Prostaglandin A₂

Item No. 10290

CAS Registry No: 130209-80-2
Formal Name: 9-oxo-15R-hydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,10-dien-1-oic acid
Synonym: 17-phenyl trinor-13,14-dihydro PGA₂
MF: C₂₃H₃₀O₄
FW: 370.5
Purity: ≥98%
UV/Vis: λ_{max}: 205 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

For long term storage, we suggest that 17-phenyl trinor-13,14-dihydro Prostaglandin A₂ (17-phenyl trinor-13,14-dihydro PGA₂) be stored as supplied at -20°C. It should be stable for at least one year.

17-phenyl trinor-13,14-dihydro 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-13,14-dihydro PGA₂ in these solvents is approximately 100 mg/ml. 17-phenyl trinor-13,14-dihydro PGA₂ is stable for at least six months in these solvents if stored at -20°C.

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-13,14-dihydro 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-13,14-dihydro PGA₂ in PBS (pH 7.2) is approximately 800 µg/ml. Store aqueous solutions of 17-phenyl trinor-13,14-dihydro PGA₂ on ice and use within 12 hours of preparation. Although the aqueous solutions of 17-phenyl trinor-13,14-dihydro PGA₂ may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Description

17-phenyl trinor-13,14-dihydro PGA₂ is a synthetic prostaglandin analog whose biological activity has not been widely reported. The PGF_{2α} analog latanoprost, which bears the same lower side chain features, has been approved as a pharmaceutical for the treatment of glaucoma.¹

Reference

1. Alm, A., Villumsen, J., Törnquist, P., *et al.* Intraocular pressure-reducing effect of PhXA41 in patients with increased eye pressure. A one-month study. *Ophthalmology* **100**, 1312-1317 (1993).

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