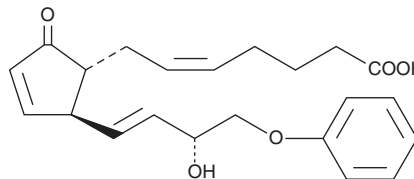


PRODUCT INFORMATION



16-phenoxy tetranor Prostaglandin A₂ Item No. 10285

CAS Registry No: 51639-10-2
Formal Name: 9-oxo-15R-hydroxy-16-phenoxy-17,18,19,20-tetranor-prosta-5Z,10,13E-trien-1-oic acid
Synonym: 16-phenoxy tetranor PGA₂
MF: C₂₂H₂₆O₅
FW: 370.4
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate
UV/Vis: λ_{max}: 219 nm



Laboratory Procedures

For long term storage, we suggest that 16-phenoxy tetranor prostaglandin A₂ (16-phenoxy tetranor PGA₂) be stored as supplied at -20°C. It should be stable for at least one year.

16-phenoxy tetranor 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 DMSO, ethanol, and dimethyl formamide purged with an inert gas can be used. The solubility of 16-phenoxy tetranor PGA₂ in these solvents is at least 100 mg/ml. 16-phenoxy tetranor 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 16-phenoxy tetranor PGA₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 16-phenoxy tetranor PGA₂ in PBS (pH 7.2) is at least 0.8 mg/ml. Store aqueous solutions of 16-phenoxy tetranor PGA₂ on ice and use within 12 hours of preparation.

Description

16-phenoxy tetranor PGA₂ is a minor metabolite found in human plasma after intravenous administration of sulprostone.¹ Its biological activity has not been studied or reported in the literature.

Reference

1. Kuhn, W., Hoyer, G., Backhus, S., *et al.* Identification of the major metabolites of the prostaglandin E₂-analogue sulprostone in human plasma, and isolation from urine (*in vivo*) and liver perfusate (*in vitro*) of female guinea-pigs. *Drug Metab. Dispos.* **19**, 920-925 (1991).

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, 08/13/2015

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