

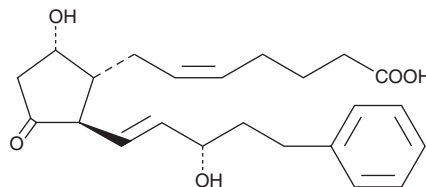
PRODUCT INFORMATION



17-phenyl trinor Prostaglandin D₂

Item No. 12810

CAS Registry No.: 85280-91-7
Formal Name: 9 α ,15S-dihydroxy-11-oxo-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-oic acid
Synonym: 17-phenyl trinor PGD₂
MF: C₂₃H₃₀O₅
FW: 386.5
Purity: \geq 98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

17-phenyl trinor Prostaglandin D₂ (17-phenyl trinor PGD₂) 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 and dimethyl formamide purged with an inert gas can be used. The solubility of 17-phenyl trinor PGD₂ in these solvents is approximately 25 and 33 mg/ml respectively. 17-phenyl trinor PGD₂ is miscible in ethanol.

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 PGD₂ 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 PGD₂ in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

17-phenyl trinor PGD₂ is a novel, chemically stable analog of PGD₂ wherein the lower side chain is modified by the addition of a phenyl group at C-17 in place of the last 3 ω -chain carbon atoms. This modification has not been reported in a D-type prostaglandin. However, in the PGE and PGF series, the analogous modification leads to a stable, somewhat more potent agonist which embodies the same biological activities as the parent prostaglandins.

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, 03/08/2024

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