

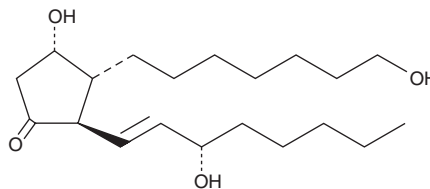
PRODUCT INFORMATION



Prostaglandin D₁ Alcohol

Item No. 12002

CAS Registry No.: 1176470-37-3
Formal Name: 1,9 α ,15S-trihydroxy-prost-13E-en-11-one
Synonym: PGD₁ Alcohol
MF: C₂₀H₃₆O₄
FW: 340.5
Purity: $\geq 99\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 1 year



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

Laboratory Procedures

Prostaglandin D₁ (PGD₁) alcohol is supplied as a crystalline solid. A stock solution may be made by dissolving the PGD₁ alcohol in the solvent of choice, which should be purged with an inert gas. PGD₁ alcohol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PGD₁ alcohol in ethanol is approximately 15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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. Organic solvent-free aqueous solutions of PGD₁ alcohol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of PGD₁ alcohol in PBS (pH 7.2) is approximately 0.075 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGD₁ alcohol is the synthetic analog of PGD₁ with a primary alcohol in place of the C-1 carboxyl. There are no known reports of its biological activity. PGD₁ is the theoretical D-series metabolite of dihomo- γ -linolenic acid (DGLA), but to date it has not been isolated as a natural product. It is an inhibitor of ADP-induced platelet aggregation in humans with an IC₅₀ value of 320 ng/ml, about 1/10 as potent as PGD₂.¹

Reference

1. Bundy, G.L., Morton, D.R., Peterson, D.C., *et al.* Synthesis and platelet aggregation inhibiting activity of prostaglandin D analogues. *J. Med. Chem.* **26(6)**, 790-799 (1983).

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, 06/06/2023

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