

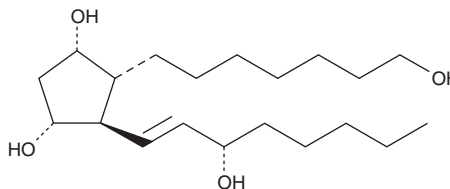
PRODUCT INFORMATION



Prostaglandin F_{1α} Alcohol

Item No. 15020

CAS Registry No.: 13487-47-3
Formal Name: 1-[3-(1-naphthalenylcarbonyl)-1H-indole-1-pentanoate]β-D-glucopyranuronic acid
Synonym: PGF_{1α} Alcohol
MF: C₂₀H₃₈O₄
FW: 342.5
Purity: ≥99%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Prostaglandin F_{1α} (PGF_{1α}) alcohol is supplied as a crystalline solid. A stock solution may be made by dissolving the PGF_{1α} alcohol in the solvent of choice, which should be purged with an inert gas. PGF_{1α} alcohol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PGF_{1α} alcohol in these solvents is approximately 10, 8, and 12 mg/ml, respectively.

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 PGF_{1α} alcohol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of PGF_{1α} alcohol in PBS (pH 7.2) is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGF_{1α} alcohol is an analog of PGF_{1α} with a primary alcohol replacing the C-1 carboxyl group. There are no published reports on the biological activity of this compound; however, the corresponding PGE₁ analog (Item No. 13020) is a relatively selective EP₃ and EP₄ receptor ligand.¹

Reference

1. Kiriya, M., Ushikubi, F., Kobayashi, T., *et al.* Ligand binding specificities of the eight types and subtypes of the mouse prostanoid receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **122**(2), 217-224 (1997).

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, 02/01/2024

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