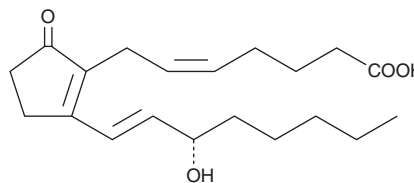


PRODUCT INFORMATION



Prostaglandin B₂ Item No. 11210

CAS Registry No.: 13367-85-6
Formal Name: 9-oxo-15S-hydroxy-prosta-5Z,8(12),13E-trien-1-oic acid
Synonym: PGB₂
MF: C₂₀H₃₀O₄
FW: 334.5
Purity: ≥98%
UV/Vis.: λ_{max}: 278 nm ε: 26,000
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

Prostaglandin B₂ (PGB₂) 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 PGB₂ in these solvents is approximately 100, 50, and 75 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. If an organic solvent-free solution of PGB₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of PGB₂ in PBS (pH 7.2) is approximately 2.0 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGB₂ is a non-enzymatic dehydration product resulting from the treatment of PGE₂ or PGA₂ with strong base. It has weak agonist activity on TP receptors and can increase pulmonary blood pressure in the rabbit at relatively high doses (5 μg/kg).¹

Reference

1. Liu, F., Orr, J.A., Wu, J. Prostaglandin B₂-induced pulmonary hypertension is mediated by TXA₂/PGH₂ receptor stimulation. *J. Am. Physiol. Soc.* **1040**, L602-L607 (1994).

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, 04/14/2020

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