

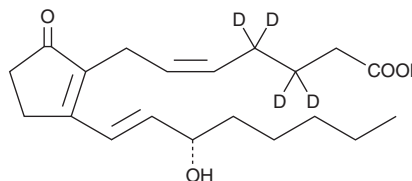
PRODUCT INFORMATION



Prostaglandin B₂-d₄

Item No. 311210

CAS Registry No.: 211105-23-6
Formal Name: 9-oxo-15S-hydroxy-prosta-5Z,8(12),13E-trien-1-oic-3,3,4,4-d₄ acid
Synonym: PGB₂-d₄
MF: C₂₀H₂₆D₄O₄
FW: 338.5
Chemical Purity: ≥99% Prostaglandin B₂
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
UV/Vis.: λ_{max}: 278 nm ε: 26,000
Supplied as: A solution in methyl acetate
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 B₂-d₄ (PGB₂-d₄) is intended for use as an internal standard for the quantification of PGB₂ (Item No. 11210) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

PGB₂-d₄ 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-d₄ in these solvents is approximately 50 mg/ml.

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., and Wu, J.-Y. 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, 04/15/2020

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