

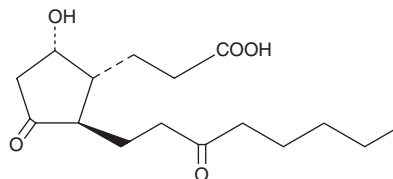
PRODUCT INFORMATION



13,14-dihydro-15-keto-tetranor Prostaglandin D₂

Item No. 13100

CAS Registry No.: 1204116-69-7
Formal Name: 9 α -hydroxy-11,15-dioxo-2,3,4,5-tetranor-prostanoic acid
Synonym: 13,14-dihydro-15-keto-tetranor PGD₂
MF: C₁₆H₂₆O₅
FW: 298.4
Purity: \geq 95%
Supplied as: A solution in ethanol
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

13,14-dihydro-15-keto-tetranor Prostaglandin D₂ (13,14-dihydro-15-keto-tetranor PGD₂) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol 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 13,14-dihydro-15-keto-tetranor PGD₂ in these solvents is approximately 50 mg/ml.

13,14-dihydro-15-keto-tetranor PGD₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 13,14-dihydro-15-keto-tetranor PGD₂ should be diluted with the aqueous buffer of choice. The solubility of 13,14-dihydro-15-keto-tetranor PGD₂ in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

A common metabolic pathway for several prostaglandins (PGs), including PGD₂, involves the reduction of the double bond between C-13 and C-14 and oxidation of the hydroxyl group at C-15, producing 13,14-dihydro-15-keto PGs. The removal of four carbons at the α -terminus and oxidation of the terminal ω -carbon produces the abundant urinary metabolites, including tetranor-PGDM.¹ 13,14-dihydro-15-keto-tetranor PGD₂ is a potential metabolite of PGD₂. It would be produced from the known metabolite 13,14-dihydro-15-keto PGD₂ (Item No. 12610), which is a known agonist for the CRTH2/DP₂ receptor.²

References

1. Song, W.L., Wang, M., Ricciotti, E., *et al.* Tetranor PGDM, an abundant urinary metabolite reflects biosynthesis of prostaglandin D₂ in mice and humans. *J. Biol. Chem.* **283**(2), 1179-1188 (2008).
2. Rangachari, P.K. and Betti, P.A. Biological activity of metabolites of PGD₂ on canine proximal colon. *Am. J. Physiol.* **264**(5 Pt 1), G886-G894 (1993).

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, 02/26/2024

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