

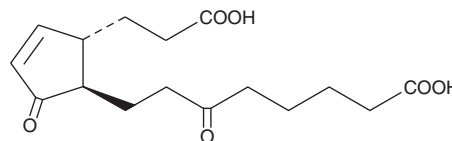
PRODUCT INFORMATION



tetranor-PGJM

Item No. 13363

CAS Registry No.: 1352751-83-7
Formal Name: 8-((1R,2S)-2-(2-carboxyethyl)-5-oxocyclopent-3-en-1-yl)-6-oxooctanoic acid
Synonym: tetranor-PGJ Metabolite
MF: C₁₆H₂₂O₆
FW: 310.3
Purity: ≥98%
UV/Vis.: λ_{max}: 219 nm
Supplied as: A solution in methyl acetate
Storage: -80°C
Stability: ≥6 months



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

Laboratory Procedures

tetranor-PGJM 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 (DMF) purged with an inert gas can be used. The solubility of tetranor-PGJM in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

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

Description

Prostaglandin D₂ (PGD₂) plays a pharmacological role in allergic and asthmatic anaphylaxis, normal physiological sleep and lowering of body temperature, as well as inhibits platelet aggregation and relaxes vascular smooth muscle.¹ tetranor-PGDM is an abundant urinary metabolite of PGD₂ that is detectable both in human and mouse and, as such, is used as a biomarker of PGD₂ biosynthesis.² tetranor-PGJM is a potential PGD₂ metabolite, formed by the elimination of the C-9 hydroxyl group. This compound may serve as a useful control in the analysis of PGD₂ biosynthesis.

References

1. Giles, H. and Leff, P. The biology and pharmacology of PGD₂. *Prostaglandins* **35**(2), 277-300 (1988).
2. 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).

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, 05/15/2023

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