

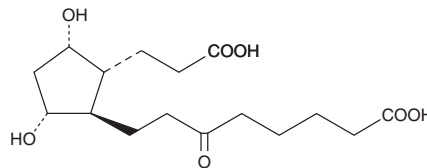
PRODUCT INFORMATION



tetranor-PGFM

Item No. 16840

CAS Registry No.: 23109-94-6
Formal Name: 9 α ,11 α -dihydroxy-15-oxo-13,14-dihydro-2,3,4,5-tetranor-prostan-1,20-dioic acid
Synonym: tetranor-Prostaglandin F Metabolite
MF: C₁₆H₂₆O₇
FW: 330.4
Purity: \geq 98%
Supplied as: A solution in methyl acetate
Storage: -80°C
Stability: \geq 1 year



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

Laboratory Procedures

tetranor-PGFM 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 tetranor-PGFM in these solvents is approximately 100 mg/ml.

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 tetranor-PGFM is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of tetranor-PGFM in PBS (pH 7.2) is approximately 10 mg/ml. Store aqueous solutions of tetranor-PGFM on ice and use within 12 hours of preparation. Although solutions of tetranor-PGFM may be stable for more than 12 hours, we recommend making a fresh preparation each day.

Description

tetranor-PGFM is the major urinary metabolite of PGF_{2 α} .¹⁻⁴ Normal healthy females excrete 7-13 μ g of tetranor-PGFM per day compared to 11-59 μ g for healthy males.² In pregnant females, tetranor-PGFM levels in the urine are 2 to 5-fold higher and diminish to pre-pregnancy levels soon after labor.²

References

1. Granström, E. and Samuelsson, B. The structure of a urinary metabolite of prostaglandin F_{2 α} in man. *J. Am. Chem. Soc.* **91**, 3398-3400 (1969).
2. Granström, E. and Samuelsson, B. On the metabolism of prostaglandin F_{2 α} in female subjects. *J. Biol. Chem.* **246**, 5254-5263 (1971).
3. Hamberg, M. Quantitative studies on prostaglandin synthesis in man III. Excretion of the major urinary metabolite of prostaglandins F_{1 α} and F_{2 α} during pregnancy. *Life Sci.* **14**, 247-252 (1974).
4. Hamberg, M. Quantitative studies on prostaglandin synthesis in man II. Determination of the major urinary metabolite of prostaglandins F_{1 α} and F_{2 α} . *Anal. Biochem.* **55**, 368-378 (1973).

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, 10/08/2018

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