

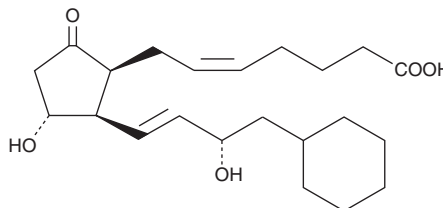
PRODUCT INFORMATION



8-iso-16-cyclohexyl-tetranor Prostaglandin E₂

Item No. 10009278

CAS Registry No.: 53319-30-5
Formal Name: 9-oxo-11 α ,15S-dihydroxy-(8 β)-16-cyclohexyl-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-oic
Synonyms: 8-iso-16-cyclohexyl-tetranor PGE₂
MF: C₂₂H₃₄O₅
FW: 378.5
Purity: \geq 97%
Supplied as: A 1 mg/ml solution in methyl acetate
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

8-iso-16-cyclohexyl-tetranor Prostaglandin E₂ (8-iso-16-cyclohexyl-tetranor PGE₂) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the 8-iso-16-cyclohexyl-tetranor PGE₂ 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 8-iso-16-cyclohexyl-tetranor PGE₂ in these solvents is approximately 50 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 8-iso-16-cyclohexyl-tetranor PGE₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 8-iso-16-cyclohexyl-tetranor PGE₂ in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

8-iso PGE₂ is one of several isoprostanes produced from polyunsaturated fatty acids during lipid peroxidation.¹ 8-iso-16-cyclohexyl-tetranor PGE₂ is a synthetic analog of 8-iso PGE₂. There are no published studies on the pharmacological properties of 8-iso-16-cyclohexyl-tetranor PGE₂.

Reference

1. Morrow, J.D., Minton, T.A., Mukundan, C.R., *et al.* Free radical-induced generation of isoprostanes *in vivo*. Evidence for the formation of D-ring and E-ring isoprostanes. *J. Biol. Chem.* **269**(6), 4317-4326 (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, 12/18/2025

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