

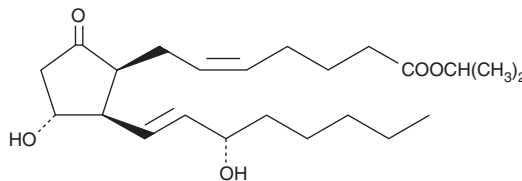
PRODUCT INFORMATION



8-iso Prostaglandin E₂ isopropyl ester

Item No. 14352

CAS Registry No.: 330589-21-4
Formal Name: (8β)-11,15S-dihydroxy-9-oxo-prosta-5Z,13E-dien-1-oic acid, 1-methylethyl ester
Synonym: 8-iso PGE₂ isopropyl ester
MF: C₂₃H₃₈O₅
FW: 394.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

8-iso Prostaglandin E₂ isopropyl ester (8-iso PGE₂ isopropyl ester) 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 8-iso PGE₂ isopropyl ester in ethanol and DMSO is approximately 50 mg/ml and approximately 30 mg/ml in DMF.

8-iso PGE₂ isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 8-iso PGE₂ isopropyl ester should be diluted with the aqueous buffer of choice. The solubility of 8-iso PGE₂ isopropyl ester in PBS (pH 7.2) is approximately 0.05 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

8-iso PGE₂ isopropyl ester is a more lipophilic form of the free acid, 8-iso PGE₂. Prostaglandin esters have enhanced lipid solubility compared to their parent compounds. They are generally hydrolyzed to the free acid upon *in vivo* administration, making the esters useful prodrugs. In general, the C-1 esters of prostaglandins show greatly diminished agonist activity *in vitro* compared to the parent free acids.

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, 02/23/2024

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