

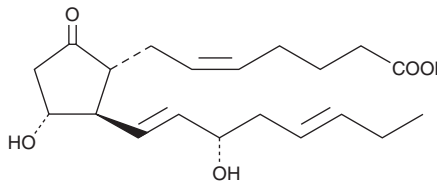
PRODUCT INFORMATION



17-*trans* Prostaglandin E₃

Item No. 14995

CAS Registry No.: 210979-33-2
Formal Name: 11 α ,15S-dihydroxy-9-oxo-prosta-5Z,13E,17E-trien-1-oic acid
Synonym: 17-*trans* PGE₃
MF: C₂₀H₃₀O₅
FW: 350.5
Purity: \geq 98%
Supplied as: A 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

17-*trans* Prostaglandin E₃ (17-*trans* PGE₃) 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 17-*trans* PGE₃ in these solvents is approximately 100 mg/ml.

17-*trans* PGE₃ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 17-*trans* PGE₃ should be diluted with the aqueous buffer of choice. The solubility of 17-*trans* PGE₃ in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

17-*trans* PGE₃ is an isomer of PGE₃ which could theoretically result from the COX metabolism of dietary trans-fatty acids.¹ There are no published reports on the biological activity of this compound.

Reference

1. Ratnayake, W.M.N., Chen, Z.Y., Pelletier, G., *et al.* Occurrence of 5c,8c,11c,15t-eicosatetraenoic acid and other unusual polyunsaturated fatty acids in rats fed partially hydrogenated canola oil. *Lipids* **29(10)**, 707-714 (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, 03/14/2024

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