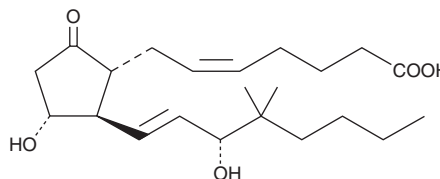


PRODUCT INFORMATION



16,16-dimethyl Prostaglandin E₂ Item No. 14750

CAS Registry No.: 39746-25-3
Formal Name: 9-oxo-11 α ,15R-dihydroxy-16,16-dimethyl-prosta-5Z,13E-dien-1-oic acid
Synonym: 16,16-dimethyl PGE₂
MF: C₂₂H₃₆O₅
FW: 380.5
Purity: \geq 95%
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

16,16-dimethyl PGE₂ 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 DMSO, ethanol, and dimethyl formamide purged with an inert gas can be used. The solubility of 16,16-dimethyl PGE₂ 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 16,16-dimethyl PGE₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 16,16-dimethyl PGE₂ in PBS (pH 7.2) is approximately 5 mg/ml. Store aqueous solutions of 16,16-dimethyl PGE₂ on ice and use within 12 hours of preparation.

Description

16,16-dimethyl PGE₂ is a competitive inhibitor of 15-hydroxy PG dehydrogenase (PGDH) but is not a substrate for the enzyme.¹ Because of its resistance to metabolism by PGDH, 16,16-dimethyl PGE₂ has a prolonged half-life *in vivo*. 16,16-dimethyl PGE₂ acts as an agonist on most EP receptor subtypes, and has been used experimentally to induce cervical ripening, uterine contraction, and prevent ulceration of the gastric mucosa in rats and dogs.^{2,3} The K_d for activation of isolated EP₂ receptors is about 1 nM.³ 16,16-dimethyl PGE₂ can be used to preserve the self-renewal properties while preventing the differentiation of hematopoietic stem cells during expansion in culture.^{4,5}

References

1. Ohno, H., Morikawa, Y., and Hirata, F. *J. Biochem.* **84**, 1485-1494 (1978).
2. Robert, A., Schultz, J.R., Nezamis, J.E., *et al. Gastroenterology* **70**, 359-370 (1976).
3. Coleman, R.A., Smith, W.L., and Narumiya, S. *Pharmacol. Rev.* **46**, 205-229 (1994).
4. Hagedorn, E.J., Durand, E.M., Fast, E.M., *et al. Exp. Cell Res.* (2014).
5. Genovese, P., Schirotti, G., Escobar, G., *et al. Nature* **510(7504)**, 235-240 (2014).

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, 12/07/2020

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