

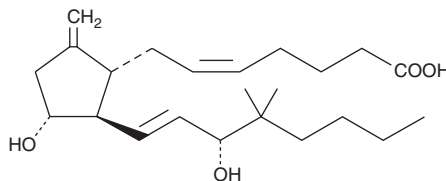
PRODUCT INFORMATION



9-deoxy-9-methylene-16,16-dimethyl Prostaglandin E₂

Item No. 14420

CAS Registry No.: 61263-35-2
Formal Name: 11 α ,15R-dihydroxy-16,16-dimethyl-9-methylene-prosta-5Z,13E-dien-1-oic acid
Synonyms: Meteneprost,
9-deoxy-9-methylene-16,16-dimethyl PGE₂
MF: C₂₃H₃₈O₄
FW: 378.6
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

9-deoxy-9-methylene-16,16-dimethyl Prostaglandin E₂ (Meteneprost) 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 meteneprost in these solvents is approximately 100 mg/ml.

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

Description

Meteneprost is a potent analog of prostaglandin E₂ (PGE₂; Item No. 14010) with an extended half life, *in vivo*. In combination with various other PG derivatives, meteneprost results in the termination of first trimester pregnancy in monkeys. A single intramuscular injection containing 0.5 mg of meteneprost and 7.5 mg of 17-phenyl trinor PGF_{1 α '} is very effective in terminating early pregnancy.¹ This PG mixture is ineffective on monkeys in their third trimester of pregnancy.¹ Meteneprost, when compared to PGE₂ and PGF_{1 α '} in monkey and rat, does not result in unwanted side effects such as fever or gastrointestinal problems.^{1,2}

References

1. Wilks, J.W. Pregnancy interception with a combination of prostaglandins: Studies in monkeys. *Science* **221(4618)**, 1407-1409 (1983).
2. Bundy, G.L., Kimball, F.A., Robert, A., *et al.* Synthesis and biological activity of 9-deoxy-9-methylene and related prostaglandins. *Adv. Prostaglandin Thromboxane Res.* **6**, 355-363 (1980).

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, 03/13/2024

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