

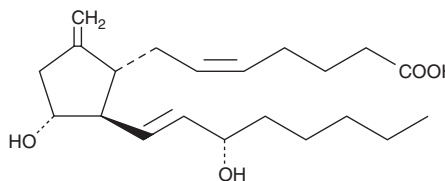
# PRODUCT INFORMATION



## 9-deoxy-9-methylene Prostaglandin E<sub>2</sub>

Item No. 14410

**CAS Registry No.:** 61263-32-9  
**Formal Name:** 11 $\alpha$ ,15S-dihydroxy-9-methylene-prosta-5Z,13E-dien-1-oic acid  
**Synonym:** 9-deoxy-9-methylene PGE<sub>2</sub>  
**MF:** C<sub>21</sub>H<sub>34</sub>O<sub>4</sub>  
**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

9-deoxy-9-methylene Prostaglandin E<sub>2</sub> (9-deoxy-9-methylene PGE<sub>2</sub>) 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 9-deoxy-9-methylene PGE<sub>2</sub> in these solvents is approximately 100 mg/ml.

9-deoxy-9-methylene PGE<sub>2</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 9-deoxy-9-methylene PGE<sub>2</sub> should be diluted with the aqueous buffer of choice. The solubility of 9-deoxy-9-methylene PGE<sub>2</sub> in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

9-deoxy-9-methylene PGE<sub>2</sub> is a synthetic analog of PGE<sub>2</sub> (Item No. 14010). 9-deoxy-9-methylene PGE<sub>2</sub> retains the biological profile of PGE<sub>2</sub> with increased stability. In the rat, 9-deoxy-9-methylene PGE<sub>2</sub> is equipotent to PGE<sub>2</sub> in decreasing blood pressure. It also stimulates the gerbil colon and primate uterus at the same potency as PGE<sub>2</sub>.<sup>1</sup>

### Reference

1. 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