

# PRODUCT INFORMATION



## 17-phenyl trinor 8-iso Prostaglandin E<sub>2</sub>

Item No. 10007931

**Formal Name:** 9-oxo-11 $\alpha$ ,15S-dihydroxy-17-phenyl-18,19,20-trinor-(8 $\beta$ )-prosta-5Z,13E-dien-1-oic acid

**MF:** C<sub>23</sub>H<sub>30</sub>O<sub>5</sub>

**FW:** 386.5

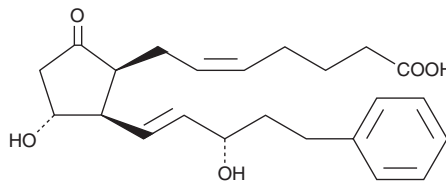
**Purity:**  $\geq$ 95%

**UV/Vis.:**  $\lambda_{\text{max}}$ : 207 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

17-phenyl trinor 8-iso Prostaglandin E<sub>2</sub> (17-phenyl trinor 8-iso PGE<sub>2</sub>) is supplied as a crystalline solid. A stock solution may be made by dissolving the 17-phenyl trinor 8-iso PGE<sub>2</sub> in the solvent of choice, which should be purged with an inert gas. 17-phenyl trinor 8-iso PGE<sub>2</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 17-phenyl trinor 8-iso PGE<sub>2</sub> in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

17-phenyl trinor 8-iso PGE<sub>2</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17-phenyl trinor 8-iso PGE<sub>2</sub> should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 17-phenyl trinor 8-iso PGE<sub>2</sub> has a solubility of approximately 0.5 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

17-phenyl trinor 8-iso PGE<sub>2</sub> is the C-8 epimer of 17-phenyl trinor PGE<sub>2</sub>, a synthetic analog of PGE<sub>2</sub>. 17-phenyl trinor PGE<sub>2</sub> is an EP<sub>1</sub> and EP<sub>3</sub> receptor agonist that causes contraction of the guinea pig ileum at a concentration of 11  $\mu$ M.<sup>1</sup> There are no published studies of the pharmacological properties of 17-phenyl trinor 8-iso PGE<sub>2</sub>.

### Reference

1. Lawrence, R.A., Jones, R.L., and Wilson, N.H. Characterization of receptors involved in the direct and indirect actions of prostaglandins E and I on the guinea-pig ileum. *Br. J. Pharmacol.* **105(2)**, 271-278 (1992).

#### 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, 04/29/2024

#### 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