

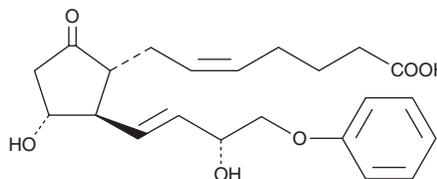
# PRODUCT INFORMATION



## 16-phenoxy tetranor Prostaglandin E<sub>2</sub>

Item No. 14760

**CAS Registry No.:** 54382-74-0  
**Formal Name:** 9-oxo-11a,15R-dihydroxy-16-phenoxy-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-oic acid  
**Synonym:** 16-phenoxy tetranor PGE<sub>2</sub>  
**MF:** C<sub>22</sub>H<sub>28</sub>O<sub>6</sub>  
**FW:** 388.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 202, 220, 271 nm  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

16-phenoxy tetranor Prostaglandin E<sub>2</sub> (16-phenoxy tetranor 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 16-phenoxy tetranor PGE<sub>2</sub> in these solvents is approximately 100 mg/ml.

16-phenoxy tetranor PGE<sub>2</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 16-phenoxy tetranor PGE<sub>2</sub> should be diluted with the aqueous buffer of choice. The solubility of 16-phenoxy tetranor PGE<sub>2</sub> in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

16-phenoxy tetranor PGE<sub>2</sub> is the free acid form of sulprostone formed by the hydrolysis of the methylsulfonamide bond.<sup>1</sup> 16-phenoxy tetranor PGE<sub>2</sub> is a minor metabolite of sulprostone found in human plasma after parenteral administration of the drug.<sup>1</sup>

### Reference

1. Kuhnz, W., Hoyer, G., Backhus, S., *et al.* Identification of the major metabolites of the prostaglandin E<sub>2</sub>-analogue sulprostone in human plasma, and isolation from urine (*in vivo*) and liver perfusate (*in vitro*) of female guinea-pigs. *Drug Metab. Dispos.* **19**(5), 920-925 (1991).

#### 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