

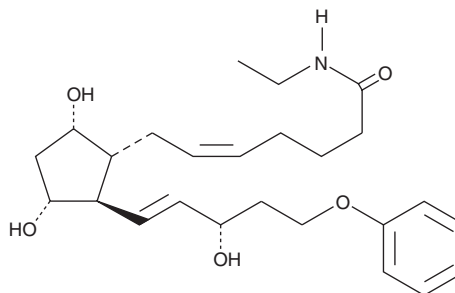
# PRODUCT INFORMATION



## 17-phenoxy trinor Prostaglandin F<sub>2α</sub> ethyl amide

Item No. 10010742

**CAS Registry No.:** 1421369-12-1  
**Formal Name:** N-ethyl-9α,11α,15S-trihydroxy-17-phenoxy-18,19,20-trinor-prosta-5Z,13E-dien-1-amide  
**Synonyms:** 17-phenoxy trinor PGF<sub>2α</sub> ethyl amide  
**MF:** C<sub>25</sub>H<sub>37</sub>NO<sub>5</sub>  
**FW:** 431.6  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

17-phenoxy trinor Prostaglandin F<sub>2α</sub> (PGF<sub>2α</sub>) ethyl amide, is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 17-phenoxy trinor PGF<sub>2α</sub> ethyl amide in these solvents is approximately 50 mg/ml.

17-phenoxy trinor PGF<sub>2α</sub> ethyl amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 17-phenoxy trinor PGF<sub>2α</sub> ethyl amide should be diluted with the aqueous buffer of choice. The solubility of 17-phenoxy trinor PGF<sub>2α</sub> ethyl amide in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PGF<sub>2α</sub>, acting through the FP receptor, causes smooth muscle contraction and exhibits potent luteolytic activity.<sup>1-3</sup> Both 17-phenyl trinor PGF<sub>2α</sub> and 16-phenoxy tetranor PGF<sub>2α</sub> are metabolically stable analogs of PGF<sub>2α</sub> and potent agonists for the FP receptor.<sup>{2058}</sup> 17-phenoxy trinor PGF<sub>2α</sub> ethyl amide is a lipophilic analog of 17-phenoxy trinor PGF<sub>2α</sub> (Item No. 10010839). Ethyl amides of PGs serve as prodrugs, as they are hydrolyzed in certain tissues to generate the bioactive free acid.

### References

1. Samuelsson, B., Goldyne, M., Granström, E., *et al.* Prostaglandins and thromboxanes. *Annu. Rev. Biochem.* **47**, 997-1029 (1978).
2. Speroff, L. and Ramwell, P.W. Prostaglandins in reproductive physiology. *Am. J. Obstet. Gynecol.* **107(7)**, 1111-1130 (1970).
3. Crankshaw, D.J. and Gaspar, V. Pharmacological characterization *in vitro* of prostanoid receptors in the myometrium of nonpregnant ewes. *J. Reprod. Fertil.* **103(1)**, 55-61 (1995).
4. Balapure, A.K., Rexroad, C.E., Jr., Kawada, K., *et al.* Structural requirements for prostaglandin analog interaction with the ovine corpus luteum prostaglandin F<sub>2α</sub> receptor. *Biochem. Pharmacol.* **38(14)**, 2375-2381 (1989).

#### 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, 04/11/2024

#### 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