

# PRODUCT INFORMATION



## 17-phenyl trinor Prostaglandin F<sub>2α</sub> cyclohexyl amide

Item No. 9000686

**Formal Name:** N-cyclohexyl-9α,11α,15S-trihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-amide

**Synonym:** Bimatoprost cyclohexyl amide

**MF:** C<sub>29</sub>H<sub>43</sub>NO<sub>4</sub>

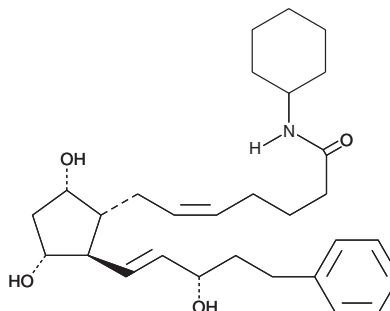
**FW:** 469.7

**Purity:** ≥95%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

17-phenyl trinor Prostaglandin F<sub>2α</sub> (17-phenyl trinor PGF<sub>2α</sub>) is supplied as a crystalline solid. A stock solution may be made by dissolving the 17-phenyl trinor PGF<sub>2α</sub> in the solvent of choice, which should be purged with an inert gas. 17-phenyl trinor PGF<sub>2α</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 17-phenyl trinor PGF<sub>2α</sub> in these solvents is approximately 30, 15, and 10 mg/ml, respectively.

17-phenyl trinor PGF<sub>2α</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17-phenyl trinor PGF<sub>2α</sub> should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 17-phenyl trinor PGF<sub>2α</sub> has a solubility of approximately 0.1 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 PGF<sub>2α</sub> is a metabolically stable, potent agonist of the PGF (FP) receptor.<sup>1,2</sup> FP agonists can be effective in modulating intraocular pressure, luteolysis, and parturition.<sup>3-5</sup> 17-phenyl trinor PGF<sub>2α</sub> cyclohexyl amide is a form of 17-phenyl trinor PGF<sub>2α</sub> with a modification of the α carbon. The addition of a cyclohexane moiety at this position strongly decreases its solubility in aqueous media, leaving it soluble solely in organic solvents. The activity of this compound at the FP receptor has not been evaluated.

### References

1. 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).
2. Lake, S., Gullberg, H., Wahlqvist, J., *et al.* Cloning of the rat and human prostaglandin F<sub>2α</sub> receptors and the expression of the rat prostaglandin F<sub>2α</sub> receptor. *FEBS Lett.* 355(3), 317-325 (1994).
3. Stjernschantz, J. and Resul, B. Phenyl substituted prostaglandin analogs for glaucoma treatment. *Drugs Future* 17(8), 691-704 (1992).
4. Dukes, M., Russell, W., and Walpole, A.L. Potent luteolytic agents related to prostaglandin F<sub>2α</sub>. *Nature* 250(464), 330-331 (1974).
5. Kobayashi, T. and Narumiya, S. Function of prostanoid receptors: Studies on knockout mice. *Prostaglandins Other Lipid Mediat.* 68-69, 557-573 (2002).

#### 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, 04/23/2024

#### 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