

PRODUCT INFORMATION



17-phenyl trinor Prostaglandin F_{2α} 1,15-lactone

Item No. 13992

Formal Name: 9α,11α,15S-trihydroxy-17-phenyl-18,19,20-trinor-prosta-5Z,13E-dien-1-oic acid, 1,15-lactone

Synonyms: Bimatoprost free acid 1,15 lactone, 17-phenyl trinor PGF_{2α} 1,15-lactone

MF: C₂₃H₃₀O₄

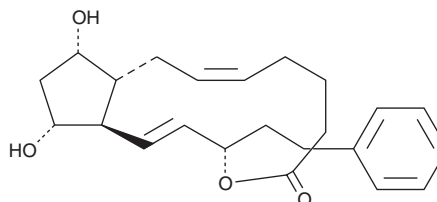
FW: 370.5

Purity: ≥98%

Supplied as: A crystalline solid

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-phenyl trinor Prostaglandin F_{2α} 1,15-lactone (17-phenyl trinor PGF_{2α} 1,15-lactone) is supplied as a crystalline solid. A stock solution may be made by dissolving the 17-phenyl trinor PGF_{2α} 1,15-lactone in the solvent of choice, which should be purged with an inert gas. 17-phenyl trinor PGF_{2α} 1,15-lactone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 17-phenyl trinor PGF_{2α} 1,15-lactone in these solvents is approximately 16, 12.5, and 14 mg/ml, respectively.

17-phenyl trinor PGF_{2α} 1,15-lactone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17-phenyl trinor PGF_{2α} 1,15-lactone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 17-phenyl trinor PGF_{2α} 1,15-lactone has a solubility of approximately 0.5 mg/ml in a 1:1 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_{2α} (Item No. 16810) is a metabolically stable analog of PGF_{2α} (Item No. 16010) with potent FP receptor agonist activity and well known intraocular pressure-reducing effects.¹⁻³ 17-phenyl trinor PGF_{2α} 1,15-lactone is the 1,15 lactone of 17-phenyl-trinor PGF_{2α}. Selective F-series PG derivatives such as this compound have been developed for suitable pharmacologic activity and an improved side-effect profile over current glaucoma therapeutics.^{4,5}

References

- Balapure, A.K., Rexroad, C.E., Jr., Kawada, K., *et al.* Structural requirements for prostaglandin analog interaction with the ovine corpus luteum prostaglandin F_{2α} receptor. *Biochem. Pharmacol.* **38(14)**, 2375-2381 (1989).
- Lake, S., Gullberg, H., Wahlqvist, J., *et al.* Cloning of the rat and human prostaglandin F_{2α} receptors and the expression of the rat prostaglandin F_{2α} receptor. *FEBS Lett.* **355(3)**, 317-325 (1994).
- Woodward, D.F., Krauss, A.H., Chen, J., *et al.* The pharmacology of bimatoprost (Lumigan™). *Surv. Ophthalmol.* **45(Suppl 4)**, S337-S345 (2001).
- Stjernschantz, J.W. From PGF_{2α}-isopropyl ester to latanoprost: A review of the development of xalatan. The proper lecture. *Invest. Ophthalmol. Vis. Sci.* **42(6)**, 1134-1145 (2001).
- Maxey, K.M. and Stanton, M.L. Internal 1,15-lactones of fluprostenol and related prostaglandin F_{2α} analogs and their use in the treatment of glaucoma and intraocular hypertension. *Cayman Chemical Company, Inc.* **US7,674,921B2** (2010).

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, 03/28/2023

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