

# PRODUCT INFORMATION



## 15-(6-nitroxyhexanoyl)-17-phenyl trinor Prostaglandin F<sub>2α</sub>

Item No. 32791

CAS Registry No.: 1194396-71-8

Formal Name: 6-(nitrooxy)-hexanoic acid, (1S,2E)-3-  
[[1R,2R,3S,5R)-2-[(2Z)-7-(ethylamino)-7-oxo-2-  
hepten-1-yl]-3,5-dihydroxycyclopentyl]-1-(2-  
phenylethyl)-2-propen-1-yl ester

Synonyms: 15-(6-nitroxyhexanoyl)-17-phenyl trinor PGF<sub>2α</sub>,  
15-(6-nitroxyhexanoyl)-Bimatoprost,  
NCX 470

MF: C<sub>31</sub>H<sub>46</sub>N<sub>2</sub>O<sub>8</sub>

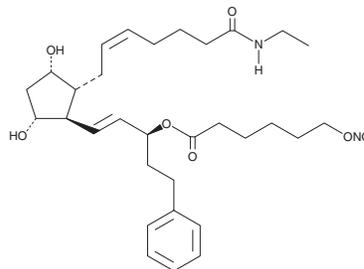
FW: 574.7

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year



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

### Laboratory Procedures

15-(6-nitroxyhexanoyl)-17-phenyl trinor Prostaglandin F<sub>2α</sub> (15-(6-nitroxyhexanoyl)-17-phenyl trinor PGF<sub>2α</sub>) 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 purged with an inert gas can be used. The solubility of 15-(6-nitroxyhexanoyl)-17-phenyl trinor PGF<sub>2α</sub> in DMSO is approximately 10 mg/ml.

### Description

15-(6-nitroxyhexanoyl)-17-phenyl trinor PGF<sub>2α</sub> is a nitric oxide-donating derivative of 17-phenyl trinor PGF<sub>2α</sub> (Item No. 16810).<sup>1</sup> It increases cGMP levels in rabbit aqueous humor and iris ciliary body when topically administered at a concentration of 0.042%. Topical administration of 15-(6-nitroxyhexanoyl)-17-phenyl trinor PGF<sub>2α</sub> (0.14%) reduces intraocular pressure (IOP) in a rabbit model of hypertonic saline-induced transient ocular hypertension. It also reduces IOP in a cynomolgus monkey model of laser-induced ocular hypertension when administered topically at a concentration of 0.042%.

### Reference

1. Impagnatiello, F., Toris, C.B., Batugo, M., *et al.* Intraocular pressure-lowering activity of NCX 470, a novel nitric oxide-donating bimatoprost in preclinical models. *Invest. Ophthalmol. Vis. Sci.* **56(11)**, 6558-6564 (2015).

#### 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, 05/06/2021

#### 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