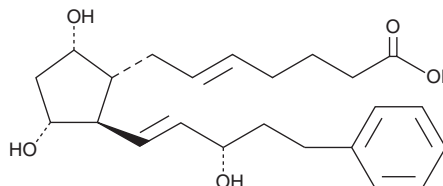


PRODUCT INFORMATION



5-*trans*-17-phenyl trinor Prostaglandin F_{2α} Item No. 9003836

CAS Registry No.: 1648894-84-1
Formal Name: (5E)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(1E,3S)-3-hydroxy-5-phenyl-1-penten-1-yl]cyclopentyl]-5-heptenoic acid
Synonyms: *trans*-Bimatoprost, 5-*trans*-17-phenyl trinor PGF_{2α}
MF: C₂₃H₃₂O₅
FW: 388.5
Purity: ≥95%
Supplied as: A 5 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

5-*trans*-17-phenyl trinor Prostaglandin F_{2α} (5-*trans*-17-phenyl trinor PGF_{2α}) is the 5-*trans* isomer of 17-phenyl trinor PGF_{2α} (Item No. 16810). It is the free acid form of 5-*trans*-17-phenyl trinor PGF_{2α} ethyl amide (Item No. 10008132) and 17-phenyl trinor PGF_{2α} dimethyl amide (Item No. 25241). 5-*trans*-17-phenyl trinor PGF_{2α} is a potential impurity in preparations of 17-phenyl trinor PGF_{2α}.¹ This item can be used as an analytical standard for the detection and quantification of 5-*trans*-17-phenyl trinor PGF_{2α} in preparations of 17-phenyl trinor PGF_{2α}.

Reference

1. Vajda, E., Hortobágyi, I., Lászlófi, I., *et al.* New process for the preparation of high purity prostaglandins. *Chinoin Gyógyszer és Vegyészeti Termékek Gyára Zrt.* **WO2015136317A1** (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, 07/14/2026

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