

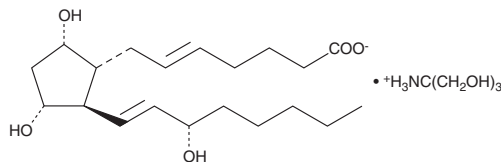
PRODUCT INFORMATION



5-*trans* Prostaglandin F_{2α} (tromethamine salt)

Item No. 16220

CAS Registry No.: 330796-56-0
Formal Name: 9α,11α,15S-trihydroxy-prosta-5E,13E-dien-1-oic acid, tris(hydroxymethyl)aminomethane salt
Synonym: 5,6-*trans* PGF_{2α} (tromethamine salt)
MF: C₂₀H₃₃O₅ • C₄H₁₂NO₃
FW: 475.6
Purity: ≥99%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that 5-*trans* Prostaglandin F_{2α} (tromethamine salt) (5-*trans* PGF_{2α} (tromethamine salt)) be stored as supplied at -20°C. It should be stable for at least two years.

5-*trans* PGF_{2α} (tromethamine salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-*trans* PGF_{2α} (tromethamine salt) in an organic solvent. 5-*trans* PGF_{2α} (tromethamine salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-*trans* PGF_{2α} (tromethamine salt) in these solvents is approximately 50 mg/ml. 5-*trans* PGF_{2α} (tromethamine salt) will be stable for at least six months in these solvents if stored at -20°C.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 5-*trans* PGF_{2α} (tromethamine salt) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 5-*trans* PGF_{2α} (tromethamine salt) in PBS is approximately 25 mg/ml compared to 10 mg/ml for 5-*trans* PGF_{2α}. We do not recommend storing the aqueous solution for more than one day.

Description

5-*trans* PGF_{2α} is the more thermodynamically stable C-5 olefin isomer of PGF_{2α} and is a common impurity in commercial lots of PGF_{2α}. There are no published reports on the biological activity of this compound.

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, 10/15/2015

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