

PRODUCT INFORMATION



6-keto Prostaglandin F_{1α} MaxSpec[®] Standard

Item No. 10007219

CAS Registry No.: 58962-34-8

Formal Name: 6-oxo-9α,11α,15S-trihydroxy-prost-13E-en-1-oic acid

Synonym: 6-keto PGF_{1α}

MF: C₂₀H₃₄O₆

FW: 370.5

Purity: ≥95%

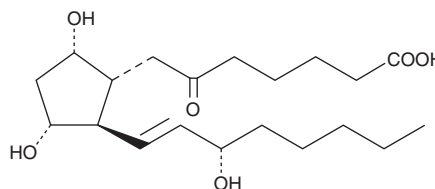
Supplied as: A solution in methyl acetate; in a deactivated glass ampule

Concentration: 100 μg/ml (nominal); see certificate of analysis for verified concentration

Storage: -20°C

Stability: ≥3 years; *Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.*

Special Conditions: Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.



Description

6-keto Prostaglandin F_{1α} (6-keto PGF_{1α}) is the inactive, non-enzymatic hydrolysis product of PGI₂ (Item No. 18220).^{1,2} 6-keto PGF_{1α} serves as a useful marker of PGI₂ biosynthesis *in vivo*.³ When [³H]-PGI₂ is injected into healthy human males, 6.6% of the radioactivity is recovered from urine as [³H]-6-keto PGF_{1α}.³

6-keto PGF_{1α} MaxSpec[®] standard is a quantitative grade standard of 6-keto PGF_{1α} (Item No. 15210) that has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. This 6-keto PGF_{1α} MaxSpec[®] standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product. **Note:** *The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.*

References

1. Pace-Asciak, C.R. Isolation, structure, and biosynthesis of 6-ketoprostaglandin F_{1α} in the rat stomach. *J. Am. Chem. Soc.* **98(8)**, 2348-2349 (1976).
2. Johnson, R.A., Morton, D.R., Kinner, J.H., et al. The chemical structure of prostaglandin X (prostacyclin). *Prostaglandins* **12(6)**, 915-928 (1976).
3. Brash, A.R., Jackson, E.K., Saggese, C.A., et al. Metabolic disposition of prostacyclin in humans. *J. Pharmacol. Exp. Ther.* **226(1)**, 78-87 (1983).

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, 09/16/2019

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