

PRODUCT INFORMATION



5 α -dihydro-11-keto Testosterone

Item No. 20200

CAS Registry No.: 32694-37-4
Formal Name: (5 α ,17 β)-17-hydroxy-androstane-3,11-dione
Synonyms: 5 α -Androstane-3,11-dione-17 β -ol, 11-keto Dihydrotestosterone, 17 β -hydroxy-5 α -Androstane-3,11-dione

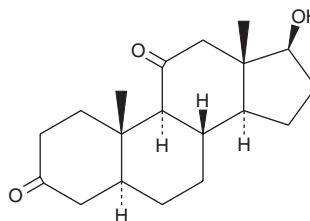
MF: C₁₉H₂₈O₃
FW: 304.4

Purity: \geq 97%

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Description

5 α -dihydro-11-keto Testosterone is an analytical reference standard that is structurally classified as a steroid hormone. It is a metabolite of 11 β -hydroxyandrostenedione and 11 β -hydroxytestosterone that has been shown to act as a full androgen receptor agonist with equal potency as that of dihydrotestosterone (Item No. 15874).¹⁻³ The androgenic activity of 5 α -dihydro-11-keto testosterone has been implicated in contributing to the androgen pool that drives castration-resistant prostate cancer cells.^{2,3} This product is intended for forensic and research applications.

References

1. Bloem, L. M., Storbeck, K. H., Schloms, L., *et al.* 11 β -Hydroxyandrostenedione returns to the steroid arena: Biosynthesis, metabolism and function. *Molecules* **18(11)**, 13228-13244 (2013).
2. Swart, A. C., and Storbeck, K. H. 11 β -Hydroxyandrostenedione: Downstream metabolism by 11 β HSD, 17 β HSD and SRD5A produces novel substrates in familiar pathway. *Mol. Cell. Endocrinol.* **408**, 114-123 (2015).
3. Storbeck, K. H., Bloem, L. M., Africander, D., *et al.* 11 β -Hydroxydihydrotestosterone and 11-ketodihydrotestosterone, novel C19 steroids with androgenic activity: A putative role in castration resistant prostate cancer? *Mol. Cell. Endocrinol.* **377(1-2)**, 135-146 (2013).

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, 11/01/2016

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