

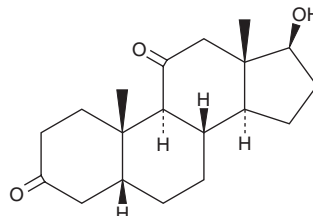
PRODUCT INFORMATION



5 β -dihydro-11-keto Testosterone

Item No. 46113

CAS Registry No.: 1420-71-9
Formal Name: (5 β)-17 β -hydroxy-androstane-3,11-dione
Synonyms: 5 β -Androstane-3,11-dione-17 β -ol, 17 β -hydroxy-5 β -Androstane-3,11-dione, 11-keto-5 β -DHT, 5 β -dihydro-11-KT
MF: C₁₉H₂₈O₃
FW: 304.4
Purity: $\geq$ 95%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

5 β -dihydro-11-keto Testosterone is supplied as a solid. A stock solution may be made by dissolving the 5 β -dihydro-11-keto testosterone in the solvent of choice, which should be purged with an inert gas. 5 β -dihydro-11-keto Testosterone is sparingly soluble (1-10 mg/ml) in methanol and chloroform.

Description

5 β -dihydro-11-keto Testosterone is a metabolite of 11-keto testosterone (Item No. 45977).¹ It is formed from 11-keto testosterone by aldo-keto reductase family 1 member D1 (AKR1D1).

Reference

1. Barnard, L., Nikolaou, N., Louw, C., *et al.* The A-ring reduction of 11-ketotestosterone is efficiently catalysed by AKR1D1 and SRD5A2 but not SRD5A1. *J. Steroid Biochem. Mol. Biol.* **202**, 105724 (2020).

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, 09/23/2026

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