

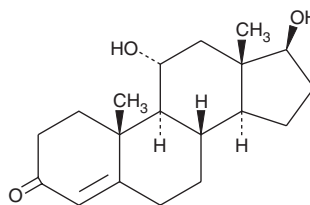
PRODUCT INFORMATION



11 α -hydroxy Testosterone

Item No. 10008647

CAS Registry No.: 3066-12-4
Formal Name: 11 α ,17 β -dihydroxy-androst-4-en-3-one
MF: C₁₉H₂₈O₃
FW: 304.4
Purity: \geq 97%
UV/Vis.: λ_{max} : 241 nm
Supplied as: A crystalline solid
Storage: 4°C
Stability: \geq 4 years



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

Laboratory Procedures

11 α -hydroxy Testosterone is supplied as a crystalline solid. A stock solution may be made by dissolving the 11 α -hydroxy testosterone in the solvent of choice, which should be purged with an inert gas. 11 α -hydroxy Testosterone is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml.

Description

11 α -hydroxy Testosterone is a metabolite of testosterone, produced by the cytochrome P450 (CYP) isoform CYP3A.¹ It can also be generated from testosterone by a mutant of CYP107A1 and from 11 α -hydroxy progesterone by the fungus *B. bassiana*.^{2,3} The biological activity of this metabolite has not been elucidated.

References

1. Zhao, Q. and Li, Z. Synthesis of 11-substituted androstenediones and testosterone as human decidual cell growth inhibitors. *Steroids* **59(3)**, 190-195 (1994).
2. Xiang, H., Tschirret-Guth, R.A., and Ortiz de Montellano, P.R. An A245T mutation conveys on cytochrome P450eryF the ability to oxidize alternative substrates. *J. Biol. Chem.* **275(46)**, 35999-36006 (2000).
3. Swizdor, A., Kolek, T., Panek, A., et al. Microbial Baeyer-Villiger oxidation of steroidal ketones using *Beauveria bassiana*: Presence of an 11 α -hydroxyl group essential to generation of D-homo lactones. *Biochim. Biophys. Acta.* **1811(4)**, 253-262 (2011).

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,10/19/2022

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