

PRODUCT INFORMATION



11-keto Testosterone-d₃

Item No. 9002754

Formal Name: 17β-hydroxy-androst-4-ene-3,11-dione-d₃

MF: C₁₉H₂₃D₃O₃

FW: 305.4

Chemical Purity: ≥98% 11-keto Testosterone

Deuterium

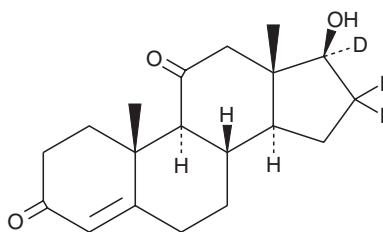
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

UV/Vis.: λ_{max}: 237 nm

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

11-keto Testosterone-d₃ (Item No. 9002754) is intended for use as an internal standard for the quantification of 11-keto testosterone (CRM) (Item No. 9002564) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

11-keto Testosterone is structurally categorized as an androgen. It is a significant hormone in certain fish, regulating the development of sexual development and behavior.^{1,2} 11-keto Testosterone binds the stickleback kidney androgen receptor (EC₅₀ = 43 nM).³ 11-keto Testosterone, at 10 nM, significantly activates the human androgen receptor expressed in zebrafish liver cells.³ This product is intended for research and forensic applications.

References

1. Yamaguchi, T., Watanuki, H., and Sakai, M. Effects of estradiol, progesterone and testosterone on the function of carp, *Cyprinus Carpio*, phagocytes *in vitro*. *Comp. Biochem. Physiol.* **129C(1)**, 49-55 (2001).
2. Godwin, J. Neuroendocrinology of sexual plasticity in teleost fishes. *Front. Neuroendocrinol.* **31(2)**, 203-216 (2010).
3. Olsson, P.-E., Berg, A.H., von Hofsten, J., *et al.* Molecular cloning and characterization of a nuclear androgen receptor activated by 11-ketotestosterone. *Reprod. Biol. Endocrinol.* **3:37**, (2005).

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, 08/16/2016

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