

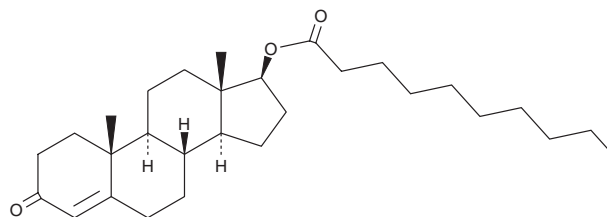
PRODUCT INFORMATION



Testosterone Decanoate

Item No. 22544

CAS Registry No.: 5721-91-5
Formal Name: 17 β -[(1-oxodecyl)oxy]-androst-4-en-3-one
Synonym: NSC 26642
MF: C₂₉H₄₆O₃
FW: 442.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 241 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Testosterone decanoate (Item No. 22544) is an analytical reference standard categorized as an anabolic androgenic steroid that can be detected in urine, blood, and hair.¹⁻³ It is an ester of the naturally occurring androgen, testosterone (Item Nos. 15645 | ISO60154) with a longer half-life. Testosterone decanoate reduces mortality and proteinuria in a murine model of lupus.⁴ Formulations containing testosterone decanoate in combination with etonogestrel (Item No. 21062) have been investigated for use as a long-acting male contraceptive.⁵ Anabolic steroids, including testosterone decanoate, have been used to enhance physical performance in racehorses and athletes.^{6,7} This product is intended for research and forensic applications.

References

1. Tretzel, L., Thomas, A., Geyer, H., *et al.* Use of dried blood spots in doping control analysis of anabolic steroid esters. *J. Pharm. Biomed. Anal.* **96**, 21-30 (2014).
2. Forsdahl, G., Erceg, D., Geisendorfer, T., *et al.* Detection of testosterone esters in blood. *Drug Test Anal.* **7(11-12)**, 983-989 (2015).
3. Kintz, P., Cirimele, V., Sachs, H., *et al.* Testing for anabolic steroids in hair from two bodybuilders. *Forensic Sci. Int.* **101(3)**, 209-216 (1999).
4. Verheul, H.A., Deckers, G.H., and Schuurs, A.H. Effects of nandrolone decanoate or testosterone decanoate on murine lupus: further evidence for a dissociation of autoimmunosuppressive and endocrine effects. *Immunopharmacology* **11(2)**, 93-99 (1986).
5. Brady, B.M., Amory, J.K., Perheentupa, A., *et al.* A multicentre study investigating subcutaneous etonogestrel implants with injectable testosterone decanoate as a potential long-acting male contraceptive. *Hum. Reprod.* **21(1)**, 284-294 (2006).
6. Forsdahl, G., Erceg, D., Geisendorfer, T., *et al.* Detection of testosterone esters in blood. *Drug Test Anal.* **7(11-12)**, 983-989 (2015).
7. You, Y., Uboh, C.E., Soma, L.R., *et al.* Simultaneous separation and determination of 16 testosterone and nandrolone esters in equine plasma using ultra high performance liquid chromatography-tandem mass spectrometry for doping control. *J. Chromatogr. A.* **1218(26)**, 3982-3993 (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, 06/15/2023

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