

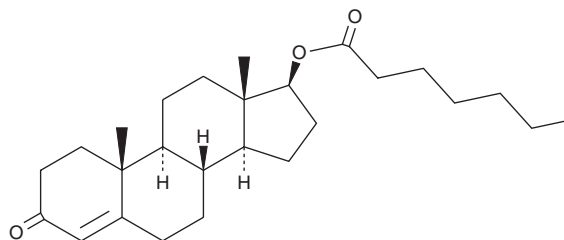
PRODUCT INFORMATION



Testosterone Enanthate

Item No. 22546

CAS Registry No.: 315-37-7
Formal Name: 17β-[(1-oxoheptyl)oxy]-androst-4-en-3-one
Synonym: NSC 17591
MF: C₂₆H₄₀O₃
FW: 400.6
Purity: ≥98%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥5 years



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

Description

Testosterone enanthate (Item No. 22546) is an analytical reference standard categorized as an anabolic androgenic steroid. It is an ester of the naturally occurring androgen, testosterone (Item Nos. 15645 | ISO60154) with a longer half-life.¹ Formulations containing testosterone enanthate have been tested for use in hypogonadism and as a potential male contraceptive.^{2,3} Formulations containing it have been associated with higher levels of total cholesterol, low density lipoproteins, and triglycerides and lower levels of high density lipoproteins.⁴ Anabolic steroids, including testosterone enanthate, have been used to enhance physical performance in racehorses and athletes, and methods to detect steroids, their derivatives, and their metabolites have been developed.^{5,6} This product is intended for research and forensic applications.

References

1. Behre, H.M. and Nieschlag, E. Testosterone Preparations for clinical use in males in *Testosterone: Action, Deficiency, Substitution*. Nieschlag, E., Behre, H.M., and Nieschlag, S., editors, 4th, Cambridge University Press, 309-335 (2012).
2. Dehennin, L. and Matsumoto, A.M. Long-term administration of testosterone enanthate to normal men: Alterations of the urinary profile of androgen metabolites potentially useful for detection of testosterone misuse in sport. *J. Steroid. Biochem. Mol. Biol.* **44(2)**, 179-189 (1993).
3. Wu, F.C.W., Balasubramanian, R., Mulders, T.M.T., et al. Oral progestogen combined with testosterone as a potential male contraceptive: Additive effects between desogestrel and testosterone enanthate in suppression of spermatogenesis, pituitary-testicular axis, and lipid metabolism. *J. Clin. Endocrinol. Metab.* **84(1)**, 112-122 (1999).
4. Jockenhövel, F., Bullmann, C., Schubert, M., et al. Influence of various modes of androgen substitution on serum lipids and lipoproteins in hypogonadal men. *Metabolism* **48(5)**, 590-596 (1999).
5. Rane, A. and Ekström, L. Androgens and doping tests: Genetic variation and pit-falls. *Br. J. Clin. Pharmacol.* **74(1)**, 3-15 (2012).
6. 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, 07/11/2024

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