

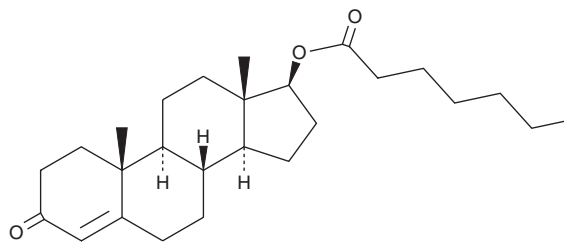
PRODUCT INFORMATION



Testosterone Enanthate (exempt preparation)

Item No. 46044

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



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

Description

Testosterone enanthate (exempt preparation) (Item No. 46044) is an analytical reference standard categorized as an anabolic androgenic steroid.^{1,2} Testosterone enanthate is a prodrug form of testosterone (Item Nos. ISO60154 | ISO00154 | 15645). It has been found in performance-enhancing drug (PED) samples seized by law enforcement.^{1,3} Formulations containing testosterone enanthate have been used in testosterone replacement therapy in patients with a condition associated with a deficiency or absence of endogenous testosterone and in the treatment of metastatic mammary cancer. Testosterone enanthate is regulated as a Schedule III compound in the United States. Testosterone enanthate (exempt preparation) (Item No. 46044) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Odoardi, S., Mestria, S., Biosa, G., *et al.* An overview on performance and image enhancing drugs (PIEDs) confiscated in Italy in the period 2017-2019. *Clin. Toxicol. (Phila)* **59(1)**, 47-52 (2021).
2. Pagonis, T.A. and Givissis, P.P. Contemporary patterns of anabolic steroids abuse and associated short term side effects in athletes. *Int. J. Orthop.* **2(4)**, 347-352 (2015).
3. Coimbra, M.A., Barbosa, F.S., Scheid, C., *et al.* Mass spectrometry determination of seized oil-based anabolic-androgenic steroids products. *Forensic Sci. Int.* **328**, 111012 (2021).

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, 05/22/2026

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