

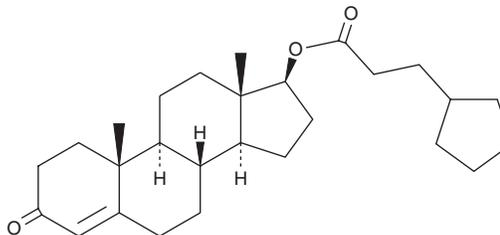
PRODUCT INFORMATION



Testosterone Cypionate

Item No. 22543

CAS Registry No.: 58-20-8
Formal Name: 17 β -(3-cyclopentyl-1-oxopropoxy)-androst-4-en-3-one
Synonym: NSC 9157
MF: C₂₇H₄₀O₃
FW: 412.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 240 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 cypionate (Item No. 22543) is an analytical reference standard categorized as an anabolic androgenic steroid. Testosterone cypionate is an ester of the naturally occurring androgen, testosterone (Item Nos. 15645 | ISO60154) with a longer half-life.¹ Formulations containing testosterone cypionate were previously used for patients with hypogonadism; however, additional esters of testosterone have been synthesized that provide improved pharmacokinetics.² Anabolic steroids, including testosterone cypionate, have been used to enhance physical performance in racehorses and athletes.^{3,4} This product is intended for research and forensic applications.

References

1. Behre, H.M., and Nieschlag, E. Testosterone preparations for clinical use in males. Testosterone: Action, deficiency, substitution (2012).
2. Behre, H.M., Abshagen, K., Oettel, M., *et al.* Intramuscular injection of testosterone undecanoate for the treatment of male hypogonadism: Phase I studies. *Eur. J. Endocrinol.* **140(5)**, 414-419 (1999).
3. Ludes, B., Kintz, P., Gromb, S., *et al.* Hair analysis of seven bodybuilders for anabolic steroids, ephedrine, and clenbuterol. *J. Forensic Sci.* **47(1)**, 211-214 (2002).
4. 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, 07/24/2023

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