

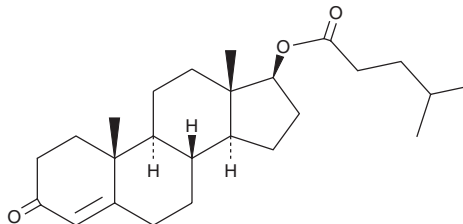
# PRODUCT INFORMATION



## Testosterone Isocaproate

Item No. 22547

**CAS Registry No.:** 15262-86-9  
**Formal Name:** (17 $\beta$ )-[(4-methyl-1-oxopentyl)oxy]-androst-4-en-3-one  
**Synonyms:** NSC 26641, Testosterone 17-Isocaproate, Testosterone 4-Methylpentanoate  
**MF:** C<sub>25</sub>H<sub>38</sub>O<sub>3</sub>  
**FW:** 386.6  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 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 isocaproate (Item No. 22547) is an analytical reference standard categorized as an anabolic androgenic steroid that can be detected in blood.<sup>1,2</sup> It is an ester of the naturally occurring androgen, testosterone (Item Nos. 15645 | ISO00154 | ISO60154) with a longer half-life.<sup>3</sup> Formulations containing testosterone isocaproate have been tested for use in hypogonadism and delayed onset puberty.<sup>4,5</sup> Anabolic steroids, including testosterone isocaproate, have been used to enhance physical performance in racehorses and athletes.<sup>6,7</sup> Testosterone isocaproate is regulated as a Schedule III compound in the United States. 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. Behre, H.M., and Nieschlag, E. Testosterone preparations for clinical use in males. Testosterone: Action, deficiency, substitution (2012).
4. Aydogdu, A., Bolu, E., Sonmez, A., *et al.* Effects of three different medications on metabolic parameters and testicular volume in patients with hypogonadotropic hypogonadism: 3-year experience. *Clin. Endocrinol. (Oxf.)* **79(2)**, 243-251 (2013).
5. Büyükgebiz, A. Treatment of constitutional delayed puberty with a combination of testosterone esters. *Horm. Res.* **44(Suppl. 3)**, 32-34 (1995).
6. 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).
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