

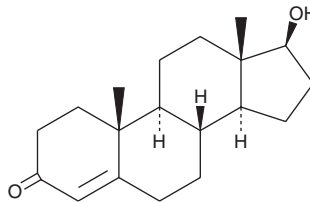
PRODUCT INFORMATION



Testosterone

Item No. 15645

CAS Registry No.: 58-22-0
Formal Name: 17 β -hydroxy-androst-4-en-3-one
Synonym: NSC 9700
MF: C₁₉H₂₈O₂
FW: 288.4
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 (Item No. 15645) is an analytical reference material categorized as an anabolic androgenic steroid.¹ Testosterone is an endogenous metabolite of androstenedione (Item No. 9001311) and estradiol (Item Nos. ISO60155 | 10006315).² Anabolic steroids, including testosterone, have been used to enhance physical performance in athletes.¹ Testosterone is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Rane, A. and Ekström, L. Androgens and doping tests: Genetic variation and pit-falls. *Br. J. Clin. Pharmacol.* **74(1)**, 3-15 (2012).
2. Baulieu, E.-E. and Mauvais-Jarvis, P. Studies on testosterone metabolism. *J. Biol. Chem.* **239(5)**, 1569-1577 (1964).

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, 06/16/2023

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