

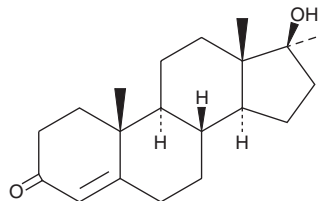
PRODUCT INFORMATION



17 α -Methyltestosterone

Item No. 24080

CAS Registry No.: 58-18-4
Formal Name: (17 β)-17-hydroxy-17-methyl-androst-4-en-3-one
Synonyms: L-589.372, NSC 139965, NSC 9701, RU-24400
MF: C₂₀H₃₀O₂
FW: 302.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 242 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

17 α -Methyltestosterone (Item No. 24080) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Anabolic androgenic steroids, including 17 α -methyltestosterone, have been used to enhance physical performance in racehorses and athletes.^{1,2} 17 α -Methyltestosterone is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Pärssinen, M. and Seppälä, T. Steroid use and long-term health risks in former athletes. *Sports Med.* **32**(2), 83-94 (2002).
2. Yu, N.H., Ho, E.N., Leung, D.K., *et al.* Screening of anabolic steroids in horse urine by liquid chromatography-tandem mass spectrometry. *J. Pharm. Biomed. Anal.* **37**(5), 1031-1038 (2005).

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, 06/21/2023

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