

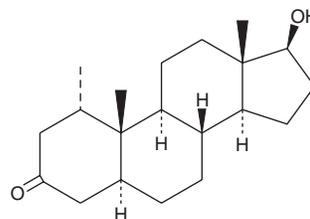
PRODUCT INFORMATION



Mesterolone

Item No. 21171

CAS Registry No.: 1424-00-6
Formal Name: (5 α)-17 β -hydroxy-1 α -methyl-androstan-3-one
Synonyms: 1 α -Methyl-5 α -androstan-17 β -ol-3-one,
1 α -methyl-5 α -Dihydrotestosterone, NSC 75054
MF: C₂₀H₃₂O₂
FW: 304.5
Purity: \geq 95%
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

Mesterolone (Item No. 21171) is an analytical reference standard categorized as an anabolic androgenic steroid. Formulations containing mesterolone have been used in patients with low testosterone or symptoms of aging male syndrome.^{1,2} Mesterolone cannot be aromatized to estrogens, unlike other androgens.^{1,3} This compound is the active ingredient in different pharmaceutical preparations that have been used medically but have also been widely applied in sports in order to improve athlete performance.⁴ Mesterolone is regulated as a Schedule III drug in the United States. This product is intended for research and forensic applications.

References

1. Carruthers, M., Cathcart, P., and Feneley, M.R. Evolution of testosterone treatment over 25 years: Symptom responses, endocrine profiles and cardiovascular changes. *Aging Male* **18(4)**, 217-227 (2015).
2. Dugeroglu, H., Ozturk, M., Atmaca, M., et al. Mesterolone treatment of aging male syndrome improves lower urinary tract symptoms. *J. Pak. Med. Assoc.* **64(12)**, 1366-1369 (2014).
3. Lowe, G.D.O., Thomson, J.E., Reavey, M.M., et al. Mesterolone: Thrombosis during treatment, and a study of its prothrombotic effects. *Br. J. Clin. Pharmacol.* **7(1)**, 107-109 (1979).
4. Dołowy, M., Pyka-Pająk, A., Filip, K., et al. A validated TLC-densitometric method for the determination of mesterolone in bulk material and in tablets. *Biomed Res. Int.* 230104 (2015).

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, 06/28/2023

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