

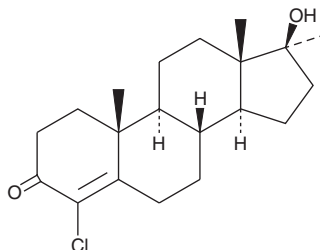
PRODUCT INFORMATION



Methylclostebol

Item No. 22448

CAS Registry No.: 5785-58-0
Formal Name: 4-chloro-17 β -hydroxy-17-methyl-androst-4-en-3-one
Synonyms: 4-chloro-17 α -Methyltestosterone, NSC 18307
MF: C₂₀H₂₉ClO₂
FW: 336.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 256 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Methylclostebol (Item No. 22448) is an analytical reference standard that is categorized as an androgenic anabolic steroid. It is a 17 α -alkylated version of clostebol (Item No. 21168; 4-chloro-17 β -hydroxy-androst-4-en-3-one), a chlorinated form of testosterone (Item Nos. ISO60154 | 15645), that can be detected in urine.^{1,2} The physiological and toxicological properties of this compound are not known. Anabolic steroids, including methylclostebol, have been used to enhance physical performance in athletes.^{2,3} Methylclostebol is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Lootens, L.M., Meuleman, P., Leroux-Roels, G., *et al.* Metabolic studies with promagnon, methylclostebol and methasterone in the uPA⁺/+SCID chimeric mice. *J. Steroid Biochem. Mol. Biol.* **127**(3-5), 374-381 (2011).
2. Franke, W.W. and Berendonk, B. Hormonal doping and androgenization of athletes: A secret program of the German democratic republic government. *Clin. Chem.* **43**(7), 1262-1279 (1997).
3. Rahnema, C.D., Crosnoe, L.E., and Kim, E.D. Designer steroids – over-the-counter supplements and their androgenic component: Review of an increasing problem. *Andrology* **3**(2), 150-155 (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, 11/11/2015

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