

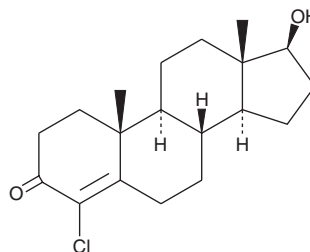
PRODUCT INFORMATION



Clotestbol

Item No. 21168

CAS Registry No.: 1093-58-9
Formal Name: 4-chloro-17 β -hydroxy-androst-4-en-3-one
Synonym: 4-chloro Testosterone
MF: C₁₉H₂₇ClO₂
FW: 322.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 254 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

Clotestbol (Item No. 21168) is an analytical reference standard that is categorized as an androgenic anabolic steroid. Clotestbol is a chlorinated form of testosterone (Item Nos. 15645 | ISO60154) that has been used as an athletic performance enhancer and to fatten cattle.¹⁻³ It is also detectible in urine. Formulations containing clotestbol have been used to treat osteoporosis, anorexia, and certain types of liver disease.⁴⁻⁶ Clotestbol is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Lu, J., Fernández-Álvarez, M., Yang, S., *et al.* New clotestbol metabolites in human urine by liquid chromatography time-of-flight tandem mass spectrometry and their application for doping control. *J. Mass. Spectrom.* **50(1)**, 191-197 (2015).
2. Debruyckere, G., de Sagher, R., and Van Peteghem, C. Clotestbol-positive urine after consumption of contaminated meat. *Clin. Chem.* **38(9)**, 1869-1873 (1992).
3. Pereira, H.M., Marques, M.A., Talhas, I.B., *et al.* Incidental clotestbol contamination in athletes after sexual intercourse. *Clin. Chem.* **50(2)**, 456-457 (2004).
4. Schanzer, W. and Donike, M. Metabolism of anabolic steroids in man: Synthesis and use of reference substances for identification of anabolic steroid metabolites. *Analytica Chim. Acta* **275(1-2)**, 23-48 (1993).
5. Calderón-Garcidueñas, L., Wen-Wang, L., Zhang, Y.J., *et al.* 8-Hydroxy-2'-deoxyguanosine, a major mutagenic oxidative DNA lesion, and DNA strand breaks in nasal respiratory epithelium of children exposed to urban pollution. *Environ. Health Perspect.* **107(6)**, 469-474 (1999).
6. Crabbe, P., Van Peteghem, C., Salden, M., *et al.* Influence of the hapten conjugation site on the characteristics of antibodies generated against metabolites of clotestbol acetate. *J. Agric. Food Chem.* **48(8)**, 3633-3638 (2000).

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, 06/22/2023

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