

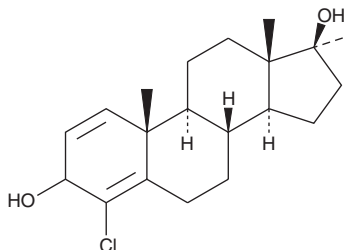
PRODUCT INFORMATION



4-Chloro-17 α -methyl-androsta-1,4-diene-3,17-diol

Item No. 22446

CAS Registry No.: 2389041-67-0
Formal Name: (17 β)-4-chloro-17-methyl-androsta-1,4-diene-3,17-diol
Synonym: Halodrol
MF: C₂₀H₂₉ClO₂
FW: 336.9
Purity: $\geq$ 95% (mixture of diastereomers)
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

4-Chloro-17 α -methyl-androsta-1,4-diene-3,17-diol (Item No. 22446) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Anabolic steroids, including 4-chloro-17 α -methyl-androsta-1,4-diene-3,17-diol, have been used to enhance physical performance in racehorses. 4-Chloro-17 α -methyl-androsta-1,4-diene-3,17-diol is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Clarke, A., Scarth, J., Teale, P., *et al.* The use of *in vitro* technologies and high-resolution/accurate-mass LC-MS to screen for metabolites of 'designer' steroids in the equine. *Drug Test Anal.* **3**(1), 74-87 (2011).

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, 11/11/205

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