

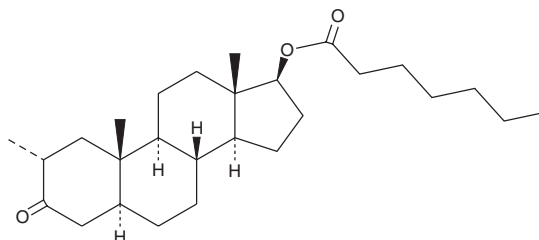
PRODUCT INFORMATION



Drostanolone Enanthate

Item No. 40656

CAS Registry No.: 13425-31-5
Formal Name: (5 α)-2 α -methyl-17 β -[(1-oxoheptyl)oxy]-androstan-3-one
Synonym: Dromostanolone Enanthate
MF: C₂₇H₄₄O₃
FW: 416.6
Purity: \geq 98%
Supplied as: A 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

Drostanolone enanthate (Item No. 40656) is an analytical reference standard categorized as an anabolic androgenic steroid prodrug of drostanolone (Item No. 21170). Drostanolone enanthate has been found in performance- and image-enhancing drug (PIED) samples seized by law enforcement.¹ Drostanolone enanthate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Odoardi, S., Mestria, S., Bioss, G., *et al.* An overview on performance and image enhancing drugs (PIEDs) confiscated in Italy in the period 2017-2019. *Clin. Toxicol. (Phila)* **59**(1), 47-52 (2021).

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, 05/28/2024

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