

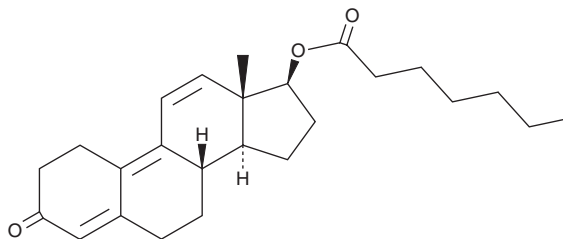
PRODUCT INFORMATION



Trenbolone Enanthate

Item No. 43900

CAS Registry No.: 1629618-98-9
Formal Name: 17 β -[(1-oxoheptyl)oxy]-estra-
4,9,11-trien-3-one
MF: C₂₅H₃₄O₃
FW: 382.5
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

Trenbolone enanthate (Item No. 43900) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Trenbolone enanthate has been found in solutions seized by government officials. Trenbolone enanthate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Coimbra, M.A., Barbosa, F.S., Scheid, C., *et al.* Mass spectrometry determination of seized oil-based anabolic-androgenic steroids products. *Forensic Sci. Int.* **328**, 111012 (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, 08/28/2025

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