

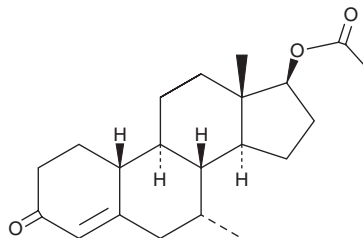
PRODUCT INFORMATION



Trestolone Acetate

Item No. 31974

CAS Registry No.: 6157-87-5
Formal Name: 17 β -(acetyloxy)-7 α -methyl-estr-4-en-3-one
Synonyms: 7 α -methyl-19-Nortestosterone Acetate, NSC 69948, U-15614
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 242 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

Trestolone acetate (Item No. 31974) is an analytical reference standard categorized as a synthetic anabolic androgenic steroid.^{1,2} Formulations containing trestolone acetate have been used to enhance physical performance in athletes.² Trestolone acetate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Noé, G., Suvisaari, J., Martin, C., *et al.* Gonadotrophin and testosterone suppression by 7 α -methyl-19-nortestosterone acetate administered by subdermal implant to healthy men. *Hum. Reprod.* **14**(9), 2200-2206 (1999).
2. Ghaphery, N.A. Performance-enhancing drugs. *Sports Med.* **26**(3), 433-442 (1995).

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, 03/27/2025

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