

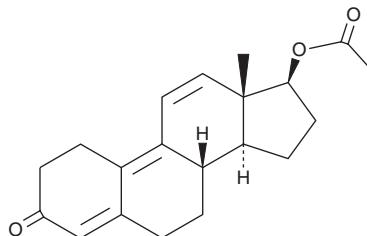
PRODUCT INFORMATION



Trenbolone Acetate

Item No. 24966

CAS Registry No.: 10161-34-9
Formal Name: 17 β -(acetyloxy)-estra-4,9,11-trien-3-one
Synonyms: RU-1697, Trenbolone 17 β -Acetate
MF: C₂₀H₂₄O₃
FW: 312.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 237, 339 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

Trenbolone acetate (Item No. 24966) is an analytical reference standard categorized as an anabolic androgenic steroid.^{1,2} Trenbolone acetate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Vernon, B.G. and Buttery, P.J. Protein turnover in rats treated with trienbolone acetate. *Br. J. Nutr.* **36(3)**, 575-579 (1976).
2. de Boer, D., Gainza Bernal, M.E., van Ooyen, R.D., *et al.* The analysis of trenbolone and the human urinary metabolites of trenbolone acetate by gas chromatography/mass spectrometry and gas chromatography/tandem mass spectrometry. *Biomol. Mass. Spectrom.* **20(8)**, 459-466 (1991).

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