

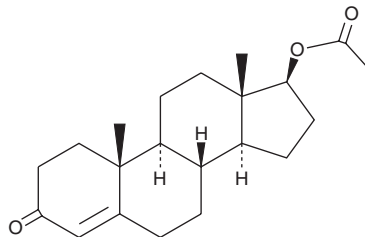
PRODUCT INFORMATION



Testosterone Acetate

Item No. 26915

CAS Registry No.: 1045-69-8
Formal Name: (17 β)-17-(acetyloxy)-androst-4-en-3-one
Synonyms: NSC 523836, 17-O-Acetyltestosterone, SKF 5647, Testosterone 17-Acetate
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 241 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

Testosterone acetate (Item No. 26915) is an analytical reference standard categorized as an anabolic androgenic steroid.^{1,2} It is an ester of the naturally occurring androgen, testosterone (Item Nos. 15645 | ISO60154). Testosterone acetate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Tretzel, L., Thomas, A., Geyer, H., *et al.* Use of dried blood spots in doping control analysis of anabolic steroid esters. *J. Pharm. Biomed. Anal.* **96**, 21-30 (2014).
2. Forsdahl, G., Erceg, D., Geisendorfer, T., *et al.* Detection of testosterone esters in blood. *Drug Test Anal.* **7(11-12)**, 983-989 (2015).

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, 07/25/2023

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