

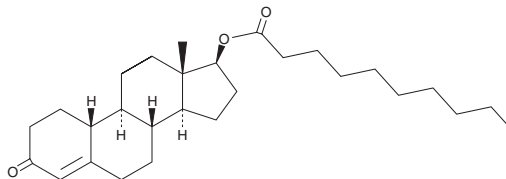
PRODUCT INFORMATION



Nandrolone Decanoate

Item No. 32546

CAS Registry No.: 360-70-3
Formal Name: 17β-[(1-oxodecyl)oxy]-estr-4-en-3-one
Synonyms: Nandrolone 17-Decanoate, Nortestosterone Decanoate
MF: C₂₈H₄₄O₃
FW: 428.7
Purity: ≥98%
UV/Vis.: λ_{max}: 240 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Nandrolone decanoate (Item No. 32546) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Nandrolone decanoate is a prodrug form of nandrolone (Item No. 21173) that has been used to enhance physical performance in athletes.² Nandrolone decanoate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Pagonis, T.A. and Givissis, P.P. Contemporary patterns of anabolic steroids abuse and associated short term side effects in athletes. *Int. J. Orthop.* **2(4)**, 347-352 (2015).
2. Berneira, L.M., de Freitas, S.C., da Silva, C.C., et al. Application of differential scanning calorimetry in the analysis of apprehended formulations of anabolic androgenic steroids. *Forensic Sci. Int.* **296**, 15-21 (2019).

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, 07/05/2023

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