

PRODUCT INFORMATION

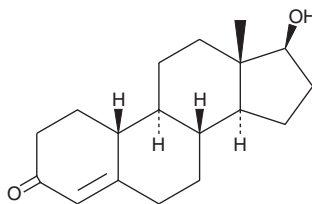


Nandrolone

Item No. 21173

CAS Registry No.: 434-22-0
Formal Name: 17 β -hydroxy-estr-4-en-3-one
Synonyms: 17 β -Nandrolone,
17 β -Nortestosterone,
19-Norandrostenolone,
19-Nortestosterone, NSC 3351

MF: C₁₈H₂₆O₂
FW: 274.4
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 207, 238, 370 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

Nandrolone (Item No. 21173) is an analytical reference standard that is structurally categorized as a steroid. It is a synthetic anabolic androgenic steroid derived from testosterone (Item No. ISO60154 | 15645). Anabolic androgenic steroids have potential therapeutic applications but are considered drugs of abuse.^{1,2} This product is intended for research and forensic applications.

References

1. Lood, Y., Eklund, A., Garle, M., *et al.* Anabolic androgenic steroids in police cases in Sweden 1999-2009. *Forensic Sci. Int.* **219**(1-3), 199-204 (2012).
2. Pan, M.M. and Kovac, J.R. Beyond testosterone cypionate: Evidence behind the use of nandrolone in male health and wellness. *Transl. Androl. Urol.* **5**(2), 213-219 (2016).

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, 06/28/2023

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