

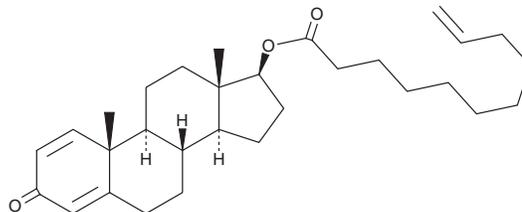
PRODUCT INFORMATION



Boldenone Undecylenate

Item No. 21632

CAS Registry No.: 13103-34-9
Formal Name: 17 β -[(1-oxo-10-undecenyl)oxy]-androsta-1,4-dien-3-one
Synonym: Boldenone 10-Undecenoate
MF: C₃₀H₄₄O₃
FW: 452.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 241 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: \geq 5 years



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

Description

Boldenone undecylenate (Item No. 21632) is an analytical reference standard that is categorized as an anabolic androgenic steroid.¹ Anabolic steroids, including boldenone undecylenate, have been used to enhance physical performance in racehorses and athletes.^{1,2} This product is intended for research and forensic applications.

References

1. Mohammed, H.H., Badawi, M.E., El-Tarabany, M.S., *et al.* Effects of boldenone undecylenate on growth performance, maintenance behaviour, reproductive hormones and carcass traits of growing rabbits. *Pol. J. Vet. Sci.* **19(2)**, 245-251 (2016).
2. O'Connor, J.J., Stillions, M.S., Reynolds, W.A., *et al.* Evaluation of boldenone undecylenate as an anabolic agent in horses. *Can. Vet. J.* **14(7)**, 154-158 (1973).

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, 07/06/2023

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