

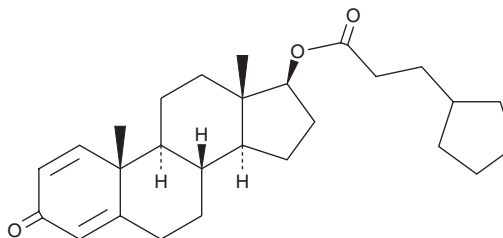
PRODUCT INFORMATION



Boldenone Cypionate

Item No. 15158

CAS Registry No.: 106505-90-2
Formal Name: 17 β -hydroxy-androsta-1,4-dien-3-one cyclopentanepropionate
MF: C₂₇H₃₈O₃
FW: 410.6
Purity: \geq 95%
UV/Vis.: λ_{max} : 244 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Description

Boldenone is an anabolic androgenic steroid and synthetic derivative of testosterone that was originally developed for veterinary use.¹ It can increase nitrogen retention, protein synthesis, and appetite, and also stimulates the release of erythropoietin in the kidneys.¹ Boldenone cypionate was synthesized as an ester of boldenone in an attempt to alter boldenone's very long half-life.² Anabolic androgenic steroid compounds such as boldenone cypionate have been used illicitly by bodybuilders and other athletes.³ This compound is intended for forensic and research purposes only.

References

1. 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).
2. 1-Dehydrotestosterone derivatives. **BE623277 19630405** (1963).
3. Socas, L., Zumbado, M., Pérez-Luzardo, O., *et al.* Hepatocellular adenomas associated with anabolic androgenic steroid abuse in bodybuilders: A report of two cases and a review of the literature. *Br. J. Sports Med.* **39(5)**, e27 (2005).

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, 04/25/2022

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