

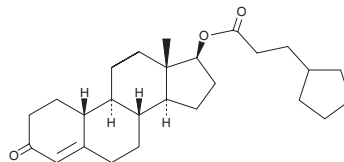
PRODUCT INFORMATION



Nandrolone Cypionate

Item No. 41907

CAS Registry No.: 601-63-8
Formal Name: 17 β -(3-cyclopentyl-1-oxopropoxy)-estr-4-en-3-one
Synonyms: NSC 40578, Nortestosterone Cypionate
MF: C₂₆H₃₈O₃
FW: 398.6
Purity: $\geq$ 98%
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: $\geq$ 4 years



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

Description

Nandrolone cypionate (Item No. 41907) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Nandrolone cypionate has been detected as an adverse analytical finding (AAF) during anti-doping testing. Nandrolone cypionate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Mullen, C., Whalley, B.J., Schifano, F., *et al.* Anabolic androgenic steroid abuse in the United Kingdom: An update. *Br. J. Pharmacol.* **177**(10), 2180-2198 (2020).

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, 09/18/2025

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