

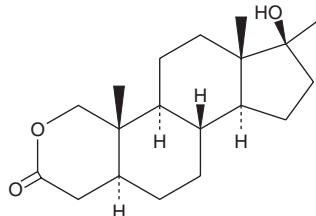
# PRODUCT INFORMATION



## Oxandrolone

Item No. 9001819

**CAS Registry No.:** 53-39-4  
**Formal Name:** (4bS,7S,9aS,9bR,11aS)-tetradecahydro-7-hydroxy-4aS,6aS,7-trimethyl-cyclopenta[5,6]naphtho[1,2-c]pyran-2(1H)-one  
**Synonyms:** NSC 67068, SC 11585, Oxandren  
**MF:** C<sub>19</sub>H<sub>30</sub>O<sub>3</sub>  
**FW:** 306.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

Oxandrolone (Item No. 9001819) is an analytical reference standard categorized as an anabolic androgenic steroid.<sup>1</sup> It has been detected as an adverse analytical finding (AAF) during anti-doping testing.<sup>2</sup> Oxandrolone is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

### References

1. Geyer, H., Schänzner, W., and Thevis, M. Anabolic agents: Recent strategies for their detection and protection from inadvertent doping. *Br. J. Sports Med.* **48(10)**, 820-826 (2014).
2. Agency, W.A.-D. 2014 Anti-doping testing figures report. (2014).

#### 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, 03/14/2019

#### 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](http://WWW.CAYMANCHEM.COM)