

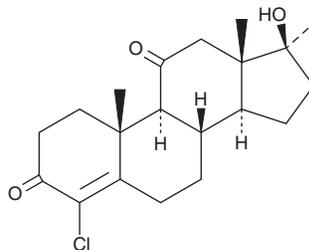
# PRODUCT INFORMATION



## 4-Chloro-17 $\beta$ -hydroxy-17-methylandro-4-ene-3,11-dione

Item No. 23310

**CAS Registry No.:** 14935-61-6  
**Formal Name:** (17 $\beta$ )-4-chloro-17-hydroxy-17-methylandro-4-ene-3,11-dione  
**MF:** C<sub>20</sub>H<sub>27</sub>ClO<sub>3</sub>  
**FW:** 350.9  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 256 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Description

4-Chloro-17 $\beta$ -hydroxy-17-methylandro-4-ene-3,11-dione (Item No. 23310) is an analytical reference standard categorized as an anabolic androgenic steroid.<sup>1</sup> It is prohibited from use in equine sports due to its anabolic enhancing properties.<sup>2</sup> 4-Chloro-17 $\beta$ -hydroxy-17-methylandro-4-ene-3,11-dione is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

### References

1. Huettenrauch, R. Synthesis of a new steroid with significant anabolic activity. *Pharmazie* **22**(3), 179 (1967).
2. Wong, A.S., Ho, E.N., Wan, T.S., *et al.* Metabolic studies of oxygundo in horses. *Anal. Chim. Acta* **891**, 190-202 (2015).

#### 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, 11/11/2015

#### 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