

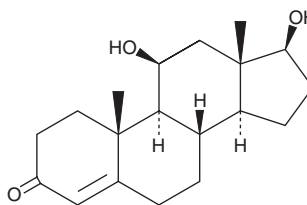
# PRODUCT INFORMATION



## 11 $\beta$ -hydroxy Testosterone

Item No. 38301

**CAS Registry No.:** 1816-85-9  
**Formal Name:** 11 $\beta$ ,17 $\beta$ -dihydroxy-androst-4-en-3-one  
**Synonym:** NSC 12168  
**MF:** C<sub>19</sub>H<sub>28</sub>O<sub>3</sub>  
**FW:** 304.4  
**Purity:**  $\geq$ 95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

11 $\beta$ -hydroxy Testosterone is supplied as a solid. A stock solution may be made by dissolving the 11 $\beta$ -hydroxy testosterone in the solvent of choice, which should be purged with an inert gas. 11 $\beta$ -hydroxy Testosterone is soluble in the organic solvent methanol.

### Description

11 $\beta$ -hydroxy Testosterone is a metabolite of testosterone.<sup>1</sup> It is formed from testosterone primarily by the cytochrome P450 (CYP) isoform CYP3A4.

### Reference

1. Choi, M.H., Skipper, P.L., Wishnok, J.S., *et al.* Characterization of testosterone 11 $\beta$ -hydroxylation catalyzed by human liver microsomal cytochromes P450. *Drug Metab. Dispos.* **33(6)**, 714-718 (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, 11/08/2023

#### 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