

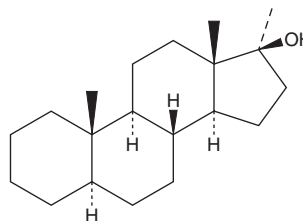
# PRODUCT INFORMATION



## 17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol

Item No. 17368

CAS Registry No.: 1229-04-5  
Formal Name: (5 $\alpha$ ,17 $\beta$ )-17-methyl-androstan-17-ol  
Synonym: Methylandrostanol  
MF: C<sub>20</sub>H<sub>34</sub>O  
FW: 290.5  
Purity:  $\geq$ 98%  
Stability:  $\geq$ 2 years at -20°C  
Supplied as: A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that 17 $\alpha$ -methyl-5 $\alpha$ -androstane-17 $\beta$ -ol be stored as supplied at -20°C. It should be stable for at least two years.

17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol is supplied as a crystalline solid. A stock solution may be made by dissolving the 17 $\alpha$ -methyl-5 $\alpha$ -androstane-17 $\beta$ -ol in the solvent of choice. 17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 17 $\alpha$ -methyl-5 $\alpha$ -androstane-17 $\beta$ -ol in ethanol is approximately 5 mg/ml and approximately 1 mg/ml in DMSO and DMF.

17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17 $\alpha$ -methyl-5 $\alpha$ -androstane-17 $\beta$ -ol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol is an analog of desoxymethyltestosterone (DMT or Madol).<sup>1</sup> 17 $\alpha$ -Methyl-5 $\alpha$ -androstane-17 $\beta$ -ol is a designer anabolic androgenic steroid that is classified as a Schedule III controlled substance in the United States. This compound is intended for forensic and research purposes only.

### Reference

1. Cawley, A.T., Blakey, K., Waller, C.C., *et al.* Detection and metabolic investigations of a novel designer steroid: 3-chloro-17 $\alpha$ -methyl-5 $\alpha$ -androstane-17 $\beta$ -ol. *Drug Test Anal.* **2015**, (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, 11/02/2015

#### 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