

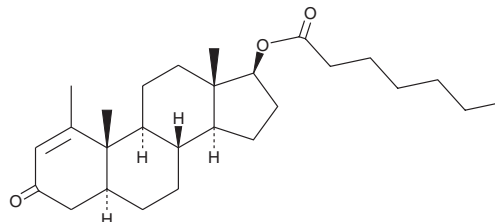
# PRODUCT INFORMATION



## Primobolan Depot

Item No. 17148

**CAS Registry No.:** 303-42-4  
**Formal Name:** (5 $\alpha$ ,17 $\beta$ )-1-methyl-17-[(1-oxoheptyl)oxy]-androst-1-en-3-one  
**Synonyms:** Methenolone Enanthate, NSC 64967, SQ 16,374  
**MF:** C<sub>27</sub>H<sub>42</sub>O<sub>3</sub>  
**FW:** 414.6  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 241 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

Primobolan depot (Item No. 17148) is an analytical reference standard that is categorized as an androgenic anabolic steroid.<sup>1</sup> Primobolan depot is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

### References

1. Nieschlag, E. and Vorona, E. Mechanisms in endocrinology: Medical consequences of doping with anabolic androgenic steroids: Effects on reproductive functions. *Eur. J. Endocrinol.* **173**(2), R47-R58 (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, 03/27/2025

#### 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