

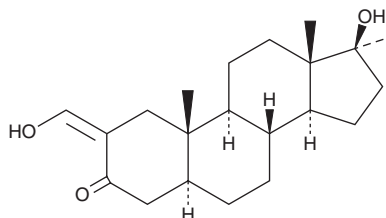
PRODUCT INFORMATION



Oxymetholone

Item No. 21643

CAS Registry No.: 434-07-1
Formal Name: (5 α ,17 β)-17-hydroxy-2-(hydroxymethylene)-17-methyl-androstan-3-one
Synonyms: C.I. 406, NSC 26198
MF: C₂₁H₃₂O₃
FW: 332.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 210, 283 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.

Laboratory Procedures

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

Oxymetholone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, oxymetholone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Oxymetholone 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

Oxymetholone is an anabolic androgenic steroid and synthetic derivative of testosterone.¹ Formulations containing oxymetholone have been used in the treatment of anemias caused by deficient red cell production.¹ Oxymetholone is regulated as a Schedule III compound in the United States.

Reference

1. Pavlatos, A.M., Fulz, O., Monberg, M.J., *et al.* Review of oxymetholone: A 17 α -alkylated anabolic-androgenic steroid. *Clin. Ther.* **23(6)**, 789-801 (2001).

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, 11/11/2025

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