

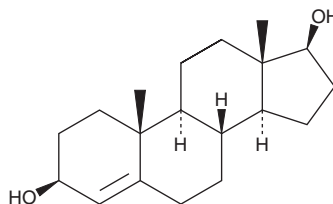
PRODUCT INFORMATION



4-Androstenediol (exempt preparation)

Item No. 46243

CAS Registry No.: 1156-92-9
Formal Name: androst-4-ene-3 β ,17 β -diol
Synonyms: 4-Adiol, 4-Androsten-3 β ,17 β -diol, NSC 12458
MF: C₁₉H₃₀O₂
FW: 290.4
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥ 3 years



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

Description

4-Androstenediol (exempt preparation) (Item No. 46243) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ 4-Androstenediol is a precursor in the biosynthesis of testosterone (Item Nos. ISO60154 | ISO00154 | 15645).² 4-Androstenediol is regulated as a Schedule III compound in the United States. 4-Androstenediol (exempt preparation) (Item No. 46243) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Poucke, C.V., Detavernier, C., Van Cauwenberghe, R., *et al.* Determination of anabolic steroids in dietary supplements by liquid chromatography-tandem mass spectrometry. *Anal. Chim. Acta* **586(1-2)**, 35-42 (2007).
2. Inaba, M. and Nakao, T. Conversion of 4-androstenediol and 5-androstenediol to testosterone, and conversion of dehydroepiandrosterone to 4-androstenediol by rat testis *in vitro*. *Endocrinol. Jpn.* **13(2)**, 160-172 (1966).

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, 06/30/2026

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