

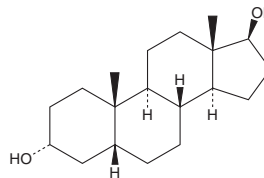
PRODUCT INFORMATION



5β-Androstane-3α,17β-diol

Item No. 33569

CAS Registry No.: 1851-23-6
Synonyms: 5β-Adiol, 5β-Androstanediol, Etiocholenediol, β-Androstane-3α,17β-diol
MF: C₁₉H₃₂O₂
FW: 292.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

5β-Androstane-3α,17β-diol is supplied as a crystalline solid. A stock solution may be made by dissolving the 5β-androstane-3α,17β-diol in the solvent of choice, which should be purged with an inert gas. 5β-Androstane-3α,17β-diol is slightly soluble in DMSO and dimethyl formamide.

Description

5β-Androstane-3α,17β-diol is a metabolite of testosterone.^{1,2} It is formed from testosterone via the 17β-hydroxyl pathway.¹

References

1. Baulieu, E.-E. and Mauvais-Jarvis, P. Studies on testosterone metabolism. *J. Biol. Chem.* **239(5)**, 1569-1577 (1964).
2. Sanyal, M.K., Orr, J.C., and Engel, L.L. The reversible reduction of testosterone to androst-4-ene-3α,17β-diol by a soluble enzyme of rat liver. *Eur. J. Biochem.* **48(1)**, 21-24 (1974).

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, 12/13/2022

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