

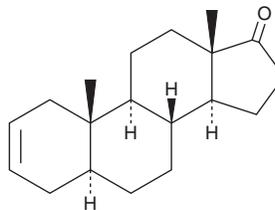
PRODUCT INFORMATION



5 α -Androst-2-ene-17-one

Item No. 21302

CAS Registry No.: 963-75-7
Formal Name: (5 α)-androst-2-en-17-one
Synonyms: NSC 44506, NSC 80614
MF: C₁₉H₂₈O
FW: 272.4
Purity: \geq 98%
Supplied as: A 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

5 α -Androst-2-ene-17-one is supplied as a solid. A stock solution may be made by dissolving the 5 α -androst-2-ene-17-one in the solvent of choice, which should be purged with an inert gas. 5 α -Androst-2-ene-17-one is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 5 α -androst-2-ene-17-one in these solvents is approximately 10 and 15 mg/ml, respectively.

Description

5 α -Androst-2-ene-17-one is a steroid and an intermediate in the synthesis of steroids and the neuromuscular blocking agent pipecuronium bromide.¹

Reference

1. Görög, S., Laukó, A., Herényi, B., et al. Analysis of steroids XXXIX. Estimation of the impurity profile of 5 α -androst-2-ene-17-one. *Chromatographia* **26**, 316-320 (1988).

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, 12/16/2022

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