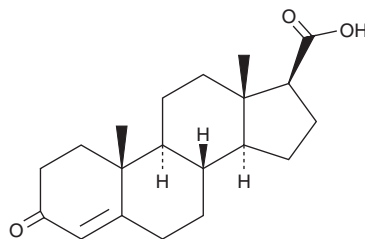


PRODUCT INFORMATION

4-Androsten-3-one-5-ene-17-carboxylic Acid

Item No. 42998

CAS Registry No.: 302-97-6
Formal Name: 3-oxo-androst-4-ene-17 β -carboxylic acid
Synonyms: AL 3793, L-552,803, NSC 226121, SKF 106224, Testosterone-17 β -carboxylic Acid
MF: C₂₀H₂₈O₃
FW: 316.4
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

4-Androsten-3-one-5-ene-17-carboxylic acid is supplied as a solid. A stock solution may be made by dissolving the 4-androsten-3-one-5-ene-17-carboxylic acid in the solvent of choice, which should be purged with an inert gas. 4-Androsten-3-one-5-ene-17-carboxylic acid is soluble ($\geq$ 10 mg/ml) in ethanol and DMSO.

Description

4-Androsten-3-one-5-ene-17-carboxylic acid is a 5 α -reductase inhibitor and a derivative of progesterone (Item No. 15876).¹ It inhibits the conversion of testosterone to dihydrotestosterone in rat ventral prostate homogenates.

Reference

1. Brooks, J.R., Baptista, E.M., Berman, C., *et al.* Response of rat ventral prostate to a new and novel 5 α -reductase inhibitor. *Endocrinology* **109**(3), 830-836 (1981).

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, 04/16/2025

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