

PRODUCT INFORMATION



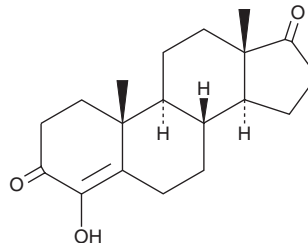
4-hydroxy Androstenedione

Item No. 18282

CAS Registry No.: 566-48-3
Formal Name: 4-hydroxy-androst-4-ene-3,17-dione
Synonym: CGP 32349, Formestane, 4-HAD, 4-Hydroxyandrostenedione, 4-Hydroxy-androst-4-ene-3,17-dione, Lentaron, NSC 282175

MF: C₁₉H₂₆O₃
FW: 302.4
Purity: ≥98%
UV/Vis.: λ_{max}: 276 nm

Storage: -20°C
Supplied as: A crystalline solid
Stability: ≥2 years



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

Laboratory Procedures

4-hydroxy Androstenedione (4-HAD) is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-HAD in the solvent of choice, which should be purged with an inert gas. 4-HAD is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 4-HAD in these solvents is approximately 10, 15, and 25 mg/ml, respectively.

4-HAD is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-HAD should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4-HAD has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-HAD is a steroidal inhibitor of aromatase (also known as cytochrome P450 19A1; K_i = 27 nM).¹ As aromatase catalyzes the conversion of androgens to estrogens, aromatase inhibitors, including 4-HAD, are used against hormone-sensitive breast cancer in menopausal women.² They are also abused in combination with anabolic steroids in racehorses and athletes.³ This product is intended for forensic and research applications.

Reference

1. Lesuisse, D., Gourvest, J.F., Hartmann, C., *et al.* Synthesis and evaluation of a new series of mechanism-based aromatase inhibitors. *J. Med. Chem.* **35(9)**, 1588-1597 (1992).
2. Smith, I.E. and Dowsett, M. Aromatase inhibitors in breast cancer. *N. Engl. J. Med.* **348(24)**, 2431-2442 (2003).
3. Cawley, A.T., Trout, G.J., Kazlauskas, R., *et al.* The detection of androstenedione abuse in sport: A mass spectrometry strategy to identify the 4-hydroxyandrostenedione metabolite. *Rapid Commun. Mass Spectrom* **22(24)**, 4147-4157 (2008).

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, 02/24/2023

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