

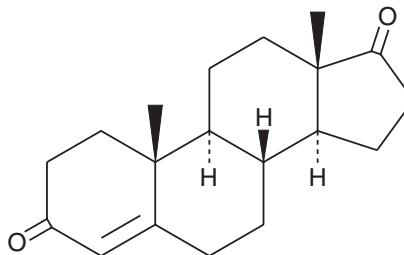
PRODUCT INFORMATION



Androstenedione

Item No. 9001311

CAS Registry No.: 63-05-8
Formal Name: androst-4-ene-3,17-dione
Synonyms: 4-Androstene-3,17-dione, Fecundin, NSC 9563, SKF 2170
MF: C₁₉H₂₆O₂
FW: 286.4
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥3 years



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

Description

Androstenedione (Item No. 9001311) is an analytical reference material categorized as an anabolic androgenic steroid.^{1,2} Androstenedione is a precursor of testosterone (Item Nos. ISO60154 | ISO00154 | 15645) and estrone (Item No. 10006485).^{2,3} Androstenedione is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Chen, F., Knecht, K., Leu, C., *et al.* Partial agonist/antagonist properties of androstenedione and 4-androsten-3 β ,17 β -diol. *J. Steroid Biochem. Mol. Biol.* **91**(4-5), 247-257 (2004).
2. Toxicology and carcinogenesis studies of androstenedione (CAS No. 63-05-8) in F344/N rats and B6C3F1 mice (gavage studies) in National Toxicology Program. National Institutes of Health, **560**, 1-171 (2010).
3. Ostojic, S.M., Calleja-Gonzalez, J., and Stojanovic, M. Steroid prohormones: Effects on body composition in athletes. *Steroids - Clinical Aspect* 11-24 (2011).

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, 02/06/2024

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