

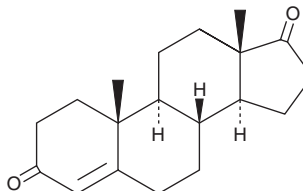
PRODUCT INFORMATION



Androstenedione (exempt preparation)

Item No. 37897

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%
UV/Vis.: λ_{max}: 243 nm
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



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

Description

Androstenedione (exempt preparation) (Item No. 37897) is an analytical reference standard categorized as an anabolic androgenic steroid.^{1,2} Androstenedione is a precursor of testosterone (Item Nos. ISO60154 | ISO00154 | 15645) and estrone (Item No. 10006485).^{1,3} Androstenedione is regulated as a Schedule III compound in the United States. Androstenedione (exempt preparation) (Item No. 37897) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Chen, F., Knecht, K., Leu, C., *et al.* Partial agonist/antagonist properties of androstenedione and 4-androsten-3β,17β-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). *Natl. Toxicol. Program Tech. Rep. Ser.* **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.* Abduljabbar, H., editor, *InTech* (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