

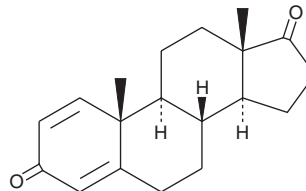
PRODUCT INFORMATION



Androsta-1,4-diene-3,17-dione

Item No. 24081

CAS Registry No.: 897-06-3
Formal Name: androsta-1,4-diene-3,17-dione
Synonyms: Boldione, NSC 49080
MF: C₁₉H₂₄O₂
FW: 284.4
Purity: ≥98%
UV/Vis.: λ_{max}: 243 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Androsta-1,4-diene-3,17-dione (Item No. 24081) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Androsta-1,4-diene-3,17-dione is a precursor in the synthesis of boldenone (Item No. 15432) and has been used to enhance athletic performance. Androsta-1,4-diene-3,17-dione is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Kim, Y., Jun, M., and Lee, W. Characterization of boldione and its metabolites in human urine by liquid chromatography/electrospray ionization mass spectrometry and gas chromatography/mass spectrometry. *Rapid Commun. Mass Spectrom.* **20(1)**, 9-20 (2006).

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, 05/20/2021

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