

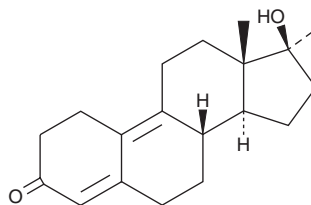
PRODUCT INFORMATION



Methyldienolone

Item No. 9002585

CAS Registry No.: 14531-89-6
Formal Name: (17β)-17-hydroxy-17-methyl-estra-4,9-dien-3-one
Synonyms: Methyl-D, Methyl-Dien, RU 3467
MF: C₁₉H₂₆O₂
FW: 286.4
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

Methyldienolone (Item No. 9002585) is an anabolic steroid that is banned by the World Anti-Doping Agency (WADA).^{1,2} It is a C-17 methylated derivative of the anabolic steroid nandrolone, which is also banned by the WADA.³ This compound is intended for forensic and research purposes only.

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

References

1. Nutting, E.F. and Calhoun, D.W. Estradienes and 2-oxaestradienes. Potent oral anabolic-androgenic agents. *Endocrinology* **84**(2), 441-442 (1969).
2. Jeong, E.S., Kim, S.-H., Cha, E.-J., *et al.* Simultaneous analysis of 210 prohibited substances in human urine by ultrafast liquid chromatography/tandem mass spectrometry in doping control. *Rapid Commun. Mass Spectrom.* **29**(4), 367-384 (2015).
3. Ayotte, C. Significance of 19-norandrosterone in athletes' urine samples. *Br. J. Sports Med.* **40 Suppl 1**, i25-i29 (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, 07/11/2019

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