

# PRODUCT INFORMATION



## Mebolazine

Item No. 9002042

**CAS Registry No.:** 3625-07-8  
**Formal Name:** [(2 $\alpha$ ,5 $\alpha$ ,17 $\beta$ )-17 $\beta$ -hydroxy-2 $\alpha$ ,17-dimethylandrostan-3-ylidene]hydrazone, 17 $\beta$ -hydroxy-2 $\alpha$ ,17-dimethyl-androstan-3-one

**Synonyms:** Dimethazine, Roxilon

**MF:** C<sub>42</sub>H<sub>68</sub>N<sub>2</sub>O<sub>2</sub>

**FW:** 633.0

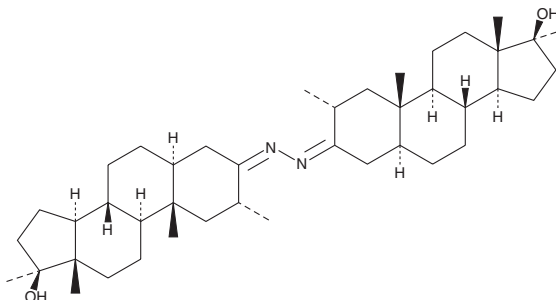
**Purity:**  $\geq$ 95%

**UV/Vis.:**  $\lambda_{\text{max}}$ : 212 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:**  $\geq$ 2 years



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

### Description

Mebolazine is a synthetic derivative of testosterone (Item No. 15645) that may bind to the androgen receptor and produce androgenic and anabolic activity.<sup>1</sup> It has been used in a variety of clinical applications to address hypoandrogenic symptoms and to generate anabolic effects; however, it carries a potential for abuse.<sup>1</sup> Mebolazine is regulated as a Schedule IV compound in the United Kingdom.<sup>2</sup> This product is intended for forensic and research applications.

### References

1. Neraud, B. and Dewailly, D. Drug-induced hyperandrogenism. *Androgen Excess Disorders in Women: Polycystic Ovary Syndrome and Other Disorders*. Azziz, R., Nestler, J.E., and Dewailly, D., editors, Humana Press, Inc. (2014).
2. Schedule 4. Part II. *Controlled drugs excepted from the prohibition on possession when in the form of a medicinal product; excluded from the application of offences arising from the prohibition on importation and exportation when imported or exported in the form of a medicinal product by any person for administration to himself; and subject to the requirements of regulations 22, 23, 26 and 27* (2014). Available from: <http://legislation.data.gov.uk/ukSI/2001/3998/schedule/4/made/data.htm?wrap=true>

#### 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, 06/17/2024

#### 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