

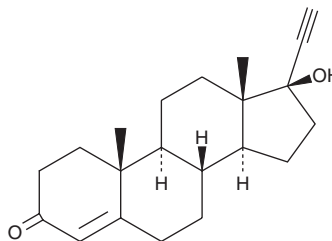
PRODUCT INFORMATION



Ethisterone

Item No. 22893

CAS Registry No.: 434-03-7
Formal Name: 17 α -hydroxy-pregn-4-en-20-yn-3-one
MF: C₂₁H₂₈O₂
FW: 312.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 241 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.

Laboratory Procedures

Ethisterone is supplied as a crystalline solid. A stock solution may be made by dissolving the ethisterone in the solvent of choice, which should be purged with an inert gas. Ethisterone is soluble in organic solvents such as ethanol, methanol, and acetonitrile. The solubility of ethisterone in these solvents is approximately 1 mg/ml.

Description

Ethisterone is an orally bioavailable synthetic progestin and a derivative of testosterone (Item Nos. ISO60154 | 15645) that binds to progesterone and androgen receptors (EC_{50} s = 23 and 23.1 nM, respectively, in yeast).¹ Ethisterone (1 μ M) downregulates progesterone receptors in MCF-7 cells.²

References

1. McRobb, L., Handelsman, D.J., Kazlauskas, R., *et al.* Structure-activity relationships of synthetic progestins in a yeast-based in vitro androgen bioassay. *J. Steroid Biochem. Mol. Biol.* **110**(1-2), 39-47 (2008).
2. Gledhill, J.M., Barker, S.A., Wanless, C., *et al.* Progesterone receptor induction by danazol in cultured cancer cells and the rat uterus. *J. Steroid. Biochem. Mol. Biol.* **43**(4), 289-296 (1992).

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, 11/12/2021

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