

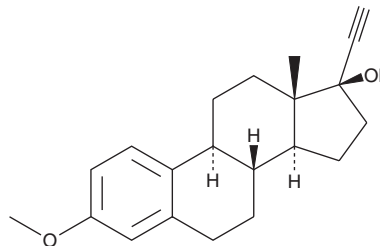
PRODUCT INFORMATION



Mestranol

Item No. 23822

CAS Registry No.: 72-33-3
Formal Name: 3-methoxy-19-norpregna-1,3,5(10)-trien-20-yn-17 α -ol
Synonyms: EE 3-ME, Ethynyl Estradiol 3-methyl ether, NSC 84032, SC-475, Δ -MVE
MF: C₂₁H₂₆O₂
FW: 310.4
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Description

Mestranol is a semisynthetic estrogen prodrug of ethynyl estradiol (Item No. 10006486), an estrogen receptor agonist.¹ It is demethylated in the liver to ethynyl estradiol. Mestranol has estrogenic properties, inducing RNA synthesis in homogenized mouse uteri following a dose of 1.6 μ g/animal, which is 16-fold higher than the dose of ethynyl estradiol needed to induce a similar response.² Formulations containing mestranol were previously used as oral contraceptives.

References

1. El-Obeid, H.A. and Al-Badr, A.A. Mestranol. *Analytical profiles of drug substances* **11**, 375-406 (1982).
2. Hahn, D.W., McGuire, J.L., Greenslade, F.C., et al. Molecular parameters involved in the estrogenicity of mestranol and ethynylestradiol. *Proc. Soc. Exp. Biol. Med.* **137(4)**, 1180-1185 (1971).

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, 11/11/2022

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