

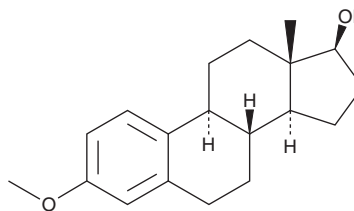
PRODUCT INFORMATION



3-O-methyl Estradiol

Item No. 42995

CAS Registry No.: 1035-77-4
Formal Name: (17 β)-3-methoxy-estra-1,3,5(10)-trien-17-ol
Synonyms: Estradiol 3-methyl ether, NSC 58851, 3-O-methyl E₂, 3-Methoxyestradiol
MF: C₁₉H₂₆O₂
FW: 286.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

3-O-methyl Estradiol is supplied as a solid. A stock solution may be made by dissolving the 3-O-methyl estradiol in the solvent of choice, which should be purged with an inert gas. 3-O-methyl Estradiol is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

3-O-methyl Estradiol is a synthetic estrogen.¹ It disrupts microtubule networks in V79 Chinese hamster cells (EC₅₀ = 9 μ M) and inhibits the proliferation of the same cells by 92% when used at a concentration of 20 μ M. 3-O-methyl Estradiol (25 mg in a 0.2 ml solution, p.o.) decreases serum cholesterol levels and increases body weight loss, indicating feminizing activity, in male rats fed a high-cholesterol diet.² It has been used as a standard for the quantification of 17 β -estradiol in hen urine.³

References

1. Aizu-Yokota, E., Susaki, A., and Sato, Y. Natural estrogens induce modulation of microtubules in Chinese hamster V79 cells in culture. *Cancer Res.* **55**(9), 1863-1868 (1995).
2. Tomarelli, R.M., Dougherty, T.M., and Bernhart, F.W. Bioassay of hypocholesterolemic steroids. *J. Pharm. Sci.* **55**(12), 1392-1395 (1966).
3. Hertelendy, F., Taylor, T.G., Mathur, R.S., *et al.* Isolation of estradiol-17 β from hen's urine and its characterization as the crystalline 3-methyl ether. *Can. J. Biochem.* **43**(9), 1379-1385 (1965).

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, 03/20/2025

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