

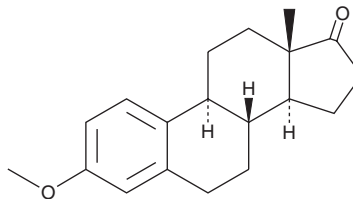
PRODUCT INFORMATION



Estrone 3-methyl ether

Item No. 27978

CAS Registry No.: 1624-62-0
Formal Name: 3-methoxy-estra-1,3,5(10)-trien-17-one
Synonyms: (+)-Estrone methyl ether,
3-O-methyl Estrone, NSC 88911
MF: C₁₉H₂₄O₂
FW: 284.4
Purity: ≥90%
Supplied as: A crystalline 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

Estrone 3-methyl ether is supplied as a crystalline solid. A stock solution may be made by dissolving the estrone 3-methyl ether in the solvent of choice, which should be purged with an inert gas. Estrone 3-methyl ether is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

Estrone 3-methyl ether is a synthetic estrogen and derivative of estrone (Item No. 10006485).¹ It stimulates proliferation of estrogen-dependent MCF-7 breast cancer cells *in vitro* when used at concentrations ranging from 0.16 to 20 μM but lacks estrogen receptor (ER) binding activity (IC₅₀s = >100 μM for ERα and ERβ). Estrone 3-methyl ether has been used in the synthesis of various estrogen receptor modulators.²

References

1. Alsayari, A., Kopel, L., Ahmed, M.S., *et al.* Design, synthesis, and biological evaluation of steroidal analogs as estrogenic/anti-estrogenic agents. *Steroids* **118**, 32-40 (2017).
2. Kuznetsov, Y.V., Levina, I.S., Scherbakov, A.M., *et al.* New estrogen receptor antagonists. 3,20-Dihydroxy-19-norpregna-1,3,5(10)-trienes: Synthesis, molecular modeling, and biological evaluation. *Eur. J. Med. Chem.* **143**, 670-682 (2018).

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, 10/05/2022

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