

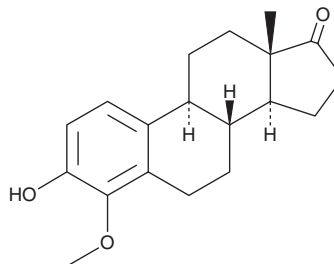
PRODUCT INFORMATION



4-methoxy Estrone

Item No. 36041

CAS Registry No.: 58562-33-7
Formal Name: 3-hydroxy-4-methoxy-estra-1,3,5(10)-trien-17-one
Synonyms: 4-MeOE₁, 4-methoxy E₁
MF: C₁₉H₂₄O₃
FW: 300.4
Purity: ≥95%
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

4-methoxy Estrone is supplied as a solid. A stock solution may be made by dissolving the 4-methoxy estrone in the solvent of choice, which should be purged with an inert gas. 4-methoxy Estrone is slightly soluble in chloroform and methanol.

Description

4-methoxy Estrone is an active metabolite of the endogenous estrogen estrone (Item No. 10006485).¹ It is formed from estrone by catechol-O-methyltransferase (COMT). 4-methoxy Estrone (10 and 100 nM) increases the proliferation of human umbilical vein endothelial cells (HUVECs).²

References

1. Dawling, S., Roodi, N., Mernaugh, R.L., *et al.* Catechol-O-methyltransferase (COMT)-mediated metabolism of catechol estrogens: Comparison of wild-type and variant COMT isoforms. *Cancer Res.* **61(18)**, 6716-6722 (2001).
2. Lippert, C., Seeger, H., Mueck, A.O., *et al.* The effects of A-ring and D-ring metabolites of estradiol on the proliferation of vascular endothelial cells. *Life Sci.* **67(13)**, 1653-1658 (2000).

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, 12/15/2022

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