

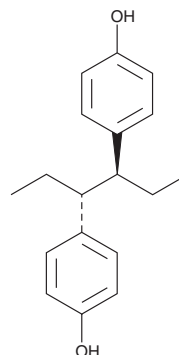
PRODUCT INFORMATION



Hexestrol

Item No. 29003

CAS Registry No.: 84-16-2
Formal Name: *rel*-4,4'-[(1*R*,2*S*)-1,2-diethyl-1,2-ethanediyl]
bis-phenol
Synonyms: Hexoestrol, *meso*-Hexestrol, NSC 9894
MF: C₁₈H₂₂O₂
FW: 270.4
Purity: ≥98%
UV/Vis.: λ_{max}: 230 nm
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

Hexestrol is supplied as a crystalline solid. A stock solution may be made by dissolving the hexestrol in the solvent of choice, which should be purged with an inert gas. Hexestrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hexestrol in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

Hexestrol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hexestrol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Hexestrol has a solubility of approximately 0.1 mg/ml in a 1:9 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hexestrol is a synthetic non-steroidal estrogen receptor ligand that binds to estrogen receptor α (ERα) and ERβ (K_ds = 0.06 and 0.06 nM for the human and rat receptors, respectively).¹ It inhibits rat liver microsomal and ox brain phospholipid liposomal lipid peroxidation (IC₅₀s = 1.5 and 2.75 μM, respectively).² It also inhibits porcine microtubule assembly and induces disassembly of preformed microtubules when used at concentrations of 50 and 100 μM, respectively, in a cell-free assay.³ Hexestrol (25 mg/animal) induces kidney tumors in male Syrian hamsters.⁴

References

1. Kuiper, G.G.J.M., Carlsson, B., Grandien, K., *et al.* Comparison of the ligand binding specificity and transcript tissue distribution of estrogen receptors α and β. *Endocrinology* **138**(3), 863-870 (1997).
2. Wiseman, H. and Halliwell, B. Carcinogenic antioxidants. Diethylstilboestrol, hexoestrol and 17 α-ethynyoestradiol. *FEBES Lett.* **332**(1-2), 159-163 (1993).
3. Sato, Y., Murai, T., Oda, T., *et al.* Inhibition of microtubule polymerization by synthetic estrogens: Formation of a ribbon structure. *J. Biochem.* **101**(5), 1247-1252 (1987).
4. Liehr, J.G., Ballatore, A.M., Dague, B.B., *et al.* Carcinogenicity and metabolic activation of hexestrol. *Chem. Biol. Interact.* **55**(1-2), 157-176 (1985).

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, 11/08/22

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