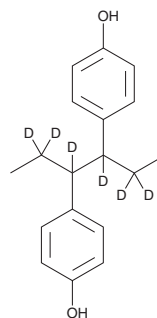


PRODUCT INFORMATION



Hexestrol-d₆ Item No. 43149

CAS Registry No.: 1219798-48-7
Formal Name: 4,4'-(hexane-3,4-diyl-2,2,3,4,5,5-d₆)diphenol
Synonyms: Hexoestrol-d₆, meso-Hexestrol-d₆
MF: C₁₈H₁₆D₆O₂
FW: 276.4
Chemical Purity: ≥98% (Hexestrol)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
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

Hexestrol-d₆ is intended for use as an internal standard for the quantification of hexestrol (Item No. 29003) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Hexestrol-d₆ is supplied as a solid. A stock solution may be made by dissolving the hexestrol-d₆ in the solvent of choice, which should be purged with an inert gas. Hexestrol-d₆ is slightly soluble (0.1-1 mg/ml) in methanol and DMSO.

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-ethynylloestradiol. *FEBS 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, 03/13/2025

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