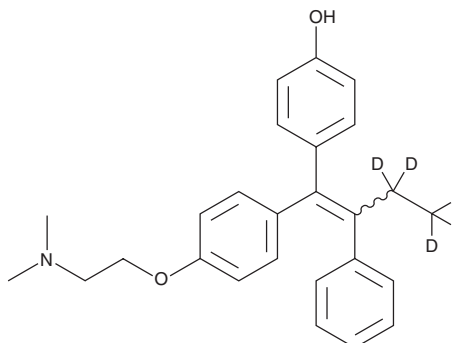


PRODUCT INFORMATION



(E/Z)-4-hydroxy Tamoxifen-d₅ Item No. 34232

CAS Registry No.: 2470232-57-4
Formal Name: 4-[1-[4-[2-(dimethylamino)ethoxy]phenyl]-2-phenyl-1-buten-1-yl]-3,3,4,4,4-d₅-phenol
Synonyms: Afimoxifene-d₅, 4-OHT-d₅
MF: C₂₆H₂₄D₅NO₂
FW: 392.6
Chemical Purity: ≥95% ((E/Z)-4-hydroxy Tamoxifen)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤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

(E/Z)-4-hydroxy Tamoxifen-d₅ is intended for use as an internal standard for the quantification of (E/Z)-4-hydroxy tamoxifen (Item No. 17308) 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).

(E/Z)-4-hydroxy Tamoxifen-d₅ is supplied as a solid. A stock solution may be made by dissolving the (E/Z)-4-hydroxy tamoxifen-d₅ in the solvent of choice, which should be purged with an inert gas. (E/Z)-4-hydroxy Tamoxifen-d₅ is soluble in organic solvents such as methanol and DMSO.

Description

(E/Z)-4-hydroxy Tamoxifen is an active metabolite of tamoxifen (Item No. 13258).¹ It is formed from tamoxifen by the cytochrome P450 (CYP) isoform CYP2D6. (E/Z)-4-hydroxy Tamoxifen is cytotoxic to MCF-7 and MDA-MB-231 breast cancer cells (IC₅₀s = 27 and 18 μM, respectively).² It also stimulates LC3 lipidation and the formation of autophagic vesicles in MCF-7 cells in a superoxide-dependent manner.³

References

1. Desta, Z., Ward, B.A., Soukhova, N.V., *et al.* Comprehensive evaluation of tamoxifen sequential biotransformation by the human cytochrome P450 system in vitro: Prominent roles for CYP3A and CYP2D6. *J. Pharmacol. Exp. Ther.* **310**(3), 1062-1075 (2004).
2. Seeger, H., Huober, J., Wallwiener, D., *et al.* Inhibition of human breast cancer cell proliferation with estradiol metabolites is as effective as with tamoxifen. *Horm. Metab. Res.* **36**(5), 277-280 (2004).
3. Duan, L., Danzer, B., Levenson, V.V., *et al.* Critical roles for nitric oxide and ERK in the completion of pro-survival autophagy in 4OHTAM-treated estrogen receptor-positive breast cancer cells. *Cancer Lett.* **353**(2), 290-300 (2014).

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, 09/08/2023

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