

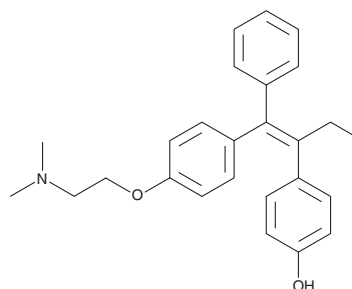
PRODUCT INFORMATION



4'-hydroxy Tamoxifen

Item No. 19519

CAS Registry No.: 82413-23-8
Formal Name: 4-[(1Z)-1-[[4-[2-(dimethylamino)ethoxy]phenyl]phenylmethylene]propyl]-phenol
MF: C₂₆H₂₉NO₂
FW: 387.5
Purity: ≥98% (may contain up to 10% of the (E) isomer)
UV/Vis.: λ_{max}: 241, 288 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

4'-hydroxy Tamoxifen is supplied as a crystalline solid. A stock solution may be made by dissolving the 4'-hydroxy tamoxifen in the solvent of choice. 4'-hydroxy Tamoxifen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 4'-hydroxy tamoxifen in ethanol and DMF is approximately 20 mg/ml and approximately 2 mg/ml in DMSO.

4'-hydroxy Tamoxifen is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4'-hydroxy tamoxifen should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 4'-hydroxy Tamoxifen has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4'-hydroxy Tamoxifen is an estrogen receptor (ER) antagonist (IC₅₀ = 2.4 nM for the rabbit receptor) and active metabolite of tamoxifen (Item Nos. 13258 | 11629).^{1,2} It is formed from tamoxifen by the cytochrome P450 (CYP) isoforms CYP2D6, CYP3A4, CYP2B6, and CYP2C19.² 4'-hydroxy Tamoxifen inhibits the proliferation of MCF-7 and MDA-MB-231 breast cancer cells (IC₅₀s = 27 and 16 μM, respectively).³ It reduces the survival percentage of HEC-1A and HEC-1B endometrial cancer cells when used at concentrations of 10 and 100 nM.⁴ Administration of liposomes encapsulating 4'-hydroxy tamoxifen (4 mg/kg twice per week) reduces tumor growth in an RPMI-8226 multiple myeloma mouse xenograft model.⁵

References

1. Löser, R., Seibel, K., Roos, W., *et al.* *In vivo* and *in vitro* antiestrogenic action of 3-hydroxytamoxifen, tamoxifen and 4-hydroxytamoxifen. *Eur. J. Cancer Clin. Oncol.* **21(8)**, 985-990 (1985).
2. Crewe, H.K., Notley, L.M., Wunsch, R.M., *et al.* Metabolism of tamoxifen by recombinant human cytochrome P450 enzymes: Formation of the 4-hydroxy, 4'-hydroxy and N-desmethyl metabolites and isomerization of trans-4-hydroxytamoxifen. *Drug Metab. Dispos.* **30(8)**, 869-874 (2002).
3. 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).
4. Cuevas, M.E. and Lindeman, T.E. *In vitro* cytotoxicity of 4'-OH-tamoxifen and estradiol in human endometrial adenocarcinoma cells HEC-1A and HEC-1B. *Oncol. Rep.* **33(1)**, 464-470 (2016).
5. Urbinati, G., Audisio, D., Marsaud, V., *et al.* Therapeutic potential of new 4-hydroxy-tamoxifen-loaded pH-gradient liposomes in a multiple myeloma experimental model. *Pharm. Res.* **27(2)**, 327-339 (2010).

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, 09/11/2023

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