

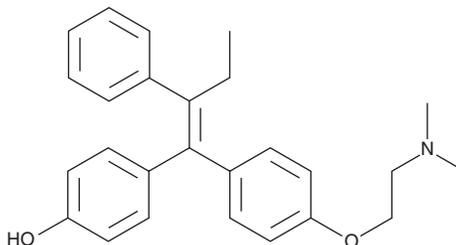
PRODUCT INFORMATION



(E)-4-hydroxy Tamoxifen

Item No. 44830

CAS Registry No.: 174592-47-3
Formal Name: 4-[(1E)-1-[4-[2-(dimethylamino)ethoxy]phenyl]-2-phenyl-1-buten-1-yl]-phenol
Synonyms: (E)-Afimoxifene, (E)-4-OHT, cis-4-hydroxy Tamoxifen
MF: C₂₆H₂₉NO₂
FW: 387.5
Purity: ≥98%
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)-4-hydroxy Tamoxifen is supplied as a solid. A stock solution may be made by dissolving the (E)-4-hydroxy tamoxifen in the solvent of choice, which should be purged with an inert gas. (E)-4-hydroxy Tamoxifen is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

(E)-4-hydroxy Tamoxifen is an active metabolite of tamoxifen (Item Nos. 13258 | 11629) and (Z)-4-hydroxy tamoxifen (Item No. 14854).¹ It is formed from tamoxifen via transisomerization of the (Z)-4-hydroxy tamoxifen intermediate by the cytochrome P450 isoforms CYP1A1, CYP1B1, CYP2B6, CYP2D6, CYP3A4, and CYP3A5. (E)-4-hydroxy Tamoxifen inhibits estradiol benzoate-induced increases in oviduct weight in quails (ED₅₀ = 2.2 mg/kg).²

Reference

1. 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).
2. Fanidi, A., Courion-Guichardaz, C., Fayard, J.M., *et al.* Effects of tamoxifen, tamoxifen metabolites, and nafoxidine on adenosine 3',5'-monophosphate phosphodiesterase: Correlations with growth inhibitory activities but not estrogen receptor affinities. *Endocrinology* **125(3)**, 1187-1193 (1989).

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, 01/23/2026

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