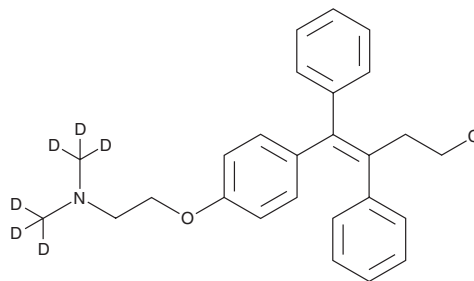


PRODUCT INFORMATION



Toremifene-d₆ Item No. 30728

Formal Name: 2-[4-[(1Z)-4-chloro-1,2-diphenyl-1-buten-1-yl]phenoxy]-N,N-di(methyl-d₃)-ethanamine
MF: C₂₆H₂₂ClD₆NO
FW: 412.0
Chemical Purity: ≥98% (Toremifene)
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

Toremifene-d₆ is intended for use as an internal standard for the quantification of toremifene (Item No. 20854) 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).

Toremifene-d₆ is supplied as a solid. A stock solution may be made by dissolving the toremifene-d₆ in the solvent of choice, which should be purged with an inert gas. Toremifene-d₆ is soluble in chloroform.

Description

Toremifene is an orally bioavailable selective estrogen receptor modulator (SERM) that evokes tissue-dependent effects similar to that of tamoxifen (Item Nos. 13258 | 11629).¹ It has been shown to block estrogen-stimulated growth of MCF-7 breast cancer cells grown in tissue culture, to inhibit ovariectomy-induced bone loss in rats, and to stimulate endometrial growth similar to that of tamoxifen.²

Reference

1. Goldstein, S.R., Siddhanti, S., Ciaccia, A.V., *et al.* A pharmacological review of selective oestrogen receptor modulators. *Hum. Reprod. Update* **6**(3), 212-224 (2000).

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, 07/20/2023

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