

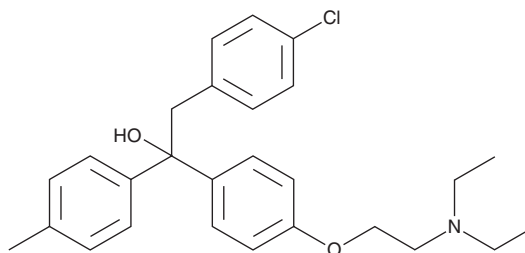
PRODUCT INFORMATION



Triparanol

Item No. 20918

CAS Registry No.: 78-41-1
Formal Name: 4-chloro- α -[4-[2-(diethylamino)ethoxy]phenyl]- α -(4-methylphenyl)-benzeneethanol
Synonyms: MER-29, NSC 65345
MF: C₂₇H₃₂ClNO₂
FW: 438.0
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 224, 274 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

Triparanol is supplied as a crystalline solid. A stock solution may be made by dissolving the triparanol in the solvent of choice. Triparanol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of triparanol in these solvents is approximately 0.125, 20, and 30 mg/ml, respectively.

Description

Triparanol is a 24-dehydro cholesterol reductase (DHCR24) inhibitor ($K_i = 0.523 \mu\text{M}$), which is an enzyme involved in the biosynthesis of cholesterol.^{1,2} It has antitumor properties, such as decreasing proliferation and inducing apoptosis in many cancer cell lines and slowing tumor growth in a mouse xenograft model.³ It can also decrease Hedgehog pathway signaling in cancer cells. Formulations containing triparanol were discontinued in the 1960s due to serious adverse side effects, including rapid cataract development.^{4,5}

References

1. Avigan, J., Steinberg, D., Vroman, H.E., *et al.* Studies of cholesterol biosynthesis. I. The identification of desmosterol in serum and tissues of animals and man treated with MER-29. *J. Biol. Chem.* **235**(11), 3123-3126 (1960).
2. Bae, S.-H., and Paik, Y.-K. Cholesterol biosynthesis from lanosterol: Development of a novel assay method and characterization of rat liver microsomal lanosterol Δ^{24} -reductase. *Biochem. J.* **326**, 609-616 (1997).
3. Bi, X., Han, X., Zhang, F., *et al.* Triparanol suppresses human tumor growth *in vitro* and *in vivo*. *Biochem. Biophys. Res. Commun.* **425**(3), 613-618 (2012).
4. Endo, A. A historical perspective on the discovery of statins. *Proc. Jpn. Acad. Ser. B. Phys. Biol. Sci.* **86**(5), 484-493 (2010).
5. Laughlin, R.C., and Carey, T.F. Cataracts in patients treated with triparanol. *JAMA* **181**(4), 339-340 (1962).

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, 02/15/2024

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