

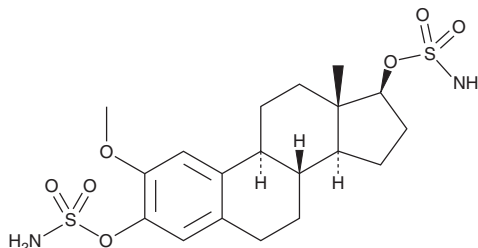
PRODUCT INFORMATION



STX140

Item No. 27863

CAS Registry No.: 401600-86-0
Formal Name: 2-methoxy-estra-1,3,5(10)-triene-3,17 β -diol, disulfamate
Synonym: 2-Methoxyestradiol-bis-sulphamate
MF: C₁₉H₂₈N₂O₇S₂
FW: 460.6
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

STX140 is supplied as a crystalline solid. A stock solution may be made by dissolving the STX140 in the solvent of choice, which should be purged with an inert gas. STX140 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of STX140 in these solvents is approximately 14, 25, and 30 mg/ml, respectively.

STX140 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, STX140 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. STX140 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

STX140 is an estrogen sulfamate with anticancer activities.¹ It inhibits steroid sulfatase with IC₅₀ values of 39 and 0.5 nM in placental microsomes and MCF-7 cancer cells, respectively. STX140 also binds to carbonic anhydrase IX and II (K_is = 70 and 270 nM, respectively).² It inhibits bovine brain tubulin assembly in a cell-free assay (IC₅₀ = 2.2 μ M) and tubule formation in human umbilical vein epithelial cells (HUVECs) when used at concentrations of 50 and 100 nM.^{3,4} STX140 inhibits proliferation of LNCaP, PC3, and MDA-MB-231 cancer cells, as well as wild-type A2780 cancer cells and adriamycin- and cisplatin-resistant A2780 cancer cells (IC₅₀s = 530, 400, 618, 330, 870, and 380 nM, respectively).^{5,6} It reduces angiogenesis in a Matrigel™ plug assay in mice and tumor growth in MCF-7 and MDA-MB-231 mouse xenograft models when used at a dose of 20 mg/kg.^{6,7}

References

1. Raobaikady, B., Purohit, A., Chander, S.K., et al. *J. Steroid Biochem. Mol. Biol.* **84**(2-3), 351-358 (2003).
2. Andring, J.T., Dohle, W., Tu, C., et al. *J. Med. Chem.* **62**(4), 2202-2212 (2019).
3. Jourdan, F., Leese, M.P., Dohle, W., et al. *J. Med. Chem.* **53**(7), 2942-2951 (2010).
4. Newman, S.P., Foster, P.A., Ho, Y.T., et al. *Br. J. Cancer* **97**(12), 1673-1682 (2007).
5. Day, J.M., Newman, S.P., Comninos, A., et al. *J. Steroid Biochem. Mol. Biol.* **84**(2-3), 317-325 (2003).
6. Foster, P.A., Ho, Y.T., Newman, S.P., et al. *Breast Cancer Res. Treat.* **111**(2), 251-260 (2008).
7. Chander, S.K., Foster, P.A., Leese, M.P., et al. *Br. J. Cancer* **96**(9), 1368-1376 (2007).

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, 06/15/2021

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