

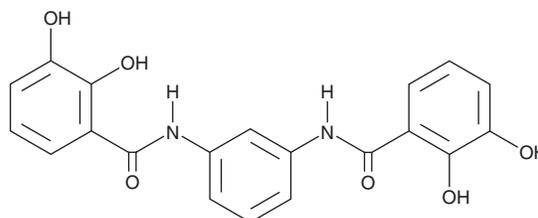
# PRODUCT INFORMATION



## MST-312

Item No. 24301

**CAS Registry No.:** 368449-04-1  
**Formal Name:** N,N'-1,3-phenylenebis[2,3-dihydroxy-benzamide]  
**Synonym:** Telomerase Inhibitor IX  
**MF:** C<sub>20</sub>H<sub>16</sub>N<sub>2</sub>O<sub>6</sub>  
**FW:** 380.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 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

MST-312 is supplied as a crystalline solid. A stock solution may be made by dissolving the MST-312 in the solvent of choice, which should be purged with an inert gas. MST-312 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MST-312 in ethanol is approximately 0.2 mg/ml and 30 mg/ml in DMSO and DMF.

MST-312 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MST-312 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MST-312 has a solubility of approximately 0.025 mg/ml in a 1:40 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MST-312 is a telomerase inhibitor (IC<sub>50</sub> = 0.67 μM in a TRAP assay).<sup>1</sup> It is selective for telomerase over DNA polymerase when used at concentrations up to 3 μM. MST-312 inhibits growth of U937 cells via telomere shortening (GI<sub>50</sub> = 1.7 μM). It reduces cell number and expression of the proliferation marker MIB-1 as well as increases DNA damage in primary pediatric ependymoma tumor cells.<sup>2</sup> MST-312 decreases survival of H460 and H1299 non-small cell lung cancer (NSCLC) cells in a dose-dependent manner with a greater effect on the aldehyde dehydrogenase positive (ALDH<sup>+</sup>) cancer stem cell population.<sup>3</sup> *In vivo*, MST-312 (40 mg/kg) reduces tumor size by 70% in an H460 mouse xenograft model. MST-312 also inhibits replication of herpes simplex virus 2 (HSV-2) and a clinical isolate of HSV-1 and decreases accumulation of early and late viral proteins in HEp-2 cells infected with HSV-1.<sup>4</sup>

### References

1. Seimiya, H., Oh-hara, T., Suzuki, T., *et al.* Telomere shortening and growth inhibition of human cancer cells by novel synthetic telomerase inhibitors MST-312, MST-295, and MST-1991. *Mol. Cancer Ther.* **1(9)**, 657-665 (2002).
2. Wong, V.C.H., Morrison, A., Tabori, U., *et al.* Telomerase inhibition as a novel therapy for pediatric ependymoma. *Brain Pathol.* **20(4)**, 780-786 (2010).
3. Serrano, D., Bleau, A.-M., Fernandez-Garcia, I., *et al.* Inhibition of telomerase activity preferentially targets aldehyde dehydrogenase-positive cancer stem-like cells in lung cancer. *Mol. Cancer* **10(96)**, (2011).
4. Haberichter, J., Roberts, S., Abbasi, I., *et al.* The telomerase inhibitor MST-312 interferes with multiple steps in the herpes simplex virus life cycle. *J. Virol.* **89(19)**, 9804-9816 (2015).

#### 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, 12/06/2022

#### 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