

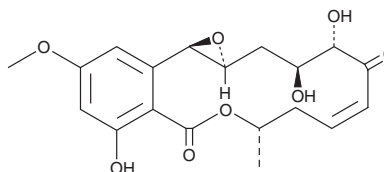
# PRODUCT INFORMATION



## Hypothemycin

Item No. 27913

**CAS Registry No.:** 76958-67-3  
**Formal Name:** (1aR,3S,4S,6Z,9S,15bR)-1a,8,9,15b-tetrahydro-3,4,12-trihydroxy-14-methoxy-9-methyl-3H-oxireno[k][2]benzoxacyclotetradecin-5,11(2H,4H)-dione  
**Synonym:** NSC 354462  
**MF:** C<sub>19</sub>H<sub>22</sub>O<sub>8</sub>  
**FW:** 378.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 221, 267, 310 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Phoma* sp.



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

### Laboratory Procedures

Hypothemycin is supplied as a solid. A stock solution may be made by dissolving the hypothemycin in the solvent of choice, which should be purged with an inert gas. Hypothemycin is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

Hypothemycin is a resorcylic acid lactone polyketide originally isolated from *H. tricothecoides* that has diverse biological activities.<sup>1-6</sup> Hypothemycin inhibits MEK (IC<sub>50</sub> = 15 nM) and other protein kinases containing a conserved cysteine residue in the ATP-binding domain, including ERK, PDGFR, VEGFR, PKD1, and MAPKAP5/MK5.<sup>3,4</sup> It also inhibits transforming growth factor β-activated kinase 1 (TAK1) *in vitro* (IC<sub>50</sub> = 33 nM).<sup>5</sup> Hypothemycin inhibits proliferation of cancer cell lines dependent on activating mutations, including A549, MV-4-11, and EOL1 cells (IC<sub>50</sub>s = 6, 0.006, and 0.0004 μM, respectively) and reduces tumor growth in Ma44 and HCT116 mouse xenograft models when administered at a dose of 25 mg/kg per day.<sup>4,6</sup>

### References

1. Agatsuma, T., Takahashi, A., Kabuto, C., *et al.* Revised structure and stereochemistry of hypothemycin. *Chem. Pharm. Bull.* **41(2)**, 373-375 (1993).
2. Xu, L., Xue, J., Wu, P., *et al.* Antifungal activity of hypothemycin against *Peronophythora litchii* in vitro and in vivo. *J. Agric. Food Chem.* **61(42)**, 10091-10095 (2013).
3. Zhao, A., Lee, S.H., Mojena, M., *et al.* Resorcylic acid lactones: Naturally occurring potent and selective inhibitors of MEK. *J. Antibiot. (Tokyo)* **52(12)**, 1086-1094 (1999).
4. Schirmer, A., Kennedy, J., Murli, S., *et al.* Targeted covalent inactivation of protein kinases by resorcylic acid lactone polyketides. *Proc. Nat. Acad. Sci. USA* **103(11)**, 4234-4239 (2006).
5. Fakhouri, L., El-Elimat, T., Hurst, D.P., *et al.* Isolation, semisynthesis, covalent docking and transforming growth factor beta-activated kinase 1 (TAK1)-inhibitory activities of (5Z)-7-oxozeaenol analogues. *Bioorg. Med. Chem.* **23(21)**, 6993-6999 (2015).
6. Tanaka, H., Nishida, K., Sugita, K., *et al.* Antitumor efficacy of hypothemycin, a new Ras-signaling inhibitor. *Jpn. J. Cancer Res.* **90(10)**, 1139-1145 (1999).

#### 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, 08/28/2023

#### 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