

# PRODUCT INFORMATION



## Cytochalasin H

Item No. 18434

**CAS Registry No.:** 53760-19-3  
**Formal Name:** (3S,3aR,4S,6S,6aR,7E,10S,12R,13E,15R,15aR)-15-(acetyloxy)-2,3,3a,4,5,6,6a,9,10,11,12,15-dodecahydro-6,12-dihydroxy-4,10,12-trimethyl-5-methylene-3-1H-cycloundec[d]isoindol-1-one

**Synonyms:** 17-Deoxo-21-acetylzygospurin D, Kodocytochalasin 1, Paspalin P1

**MF:** C<sub>30</sub>H<sub>39</sub>NO<sub>5</sub>

**FW:** 493.6

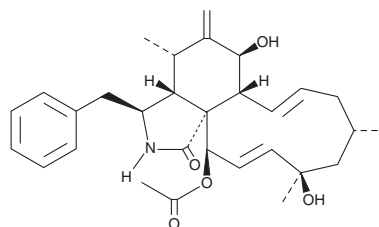
**Purity:** ≥95%

**Supplied as:** A lyophilisate

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/*Phomopsis* sp.



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

### Laboratory Procedures

Cytochalasin H is supplied as a lyophilisate. A stock solution may be made by dissolving the cytochalasin H in the solvent of choice, which should be purged with an inert gas. Cytochalasin H is soluble in organic solvents such as ethanol, methanol, DMSO and dimethyl formamide. We do not recommend storing the aqueous solution for more than one day.

### Description

The cytochalasins are cell-permeable fungal metabolites that inhibit actin polymerization.<sup>1-4</sup> This interferes with such diverse processes as cell movement, growth, phagocytosis, degranulation, and secretion.<sup>5-8</sup> Cytochalasin H is a potent inhibitor of actin incorporation into filaments.<sup>9,10</sup> Like cytochalasin D (Item No. 11330), it also increases the steady-state diffusion coefficients of filaments, suggesting that it stimulates filament shortening.<sup>9</sup> Cytochalasin H suppresses the production of reactive oxygen species by stimulated human neutrophils, blocks endocytosis of CD59 in lymphocytes, and shows anti-angiogenic activity both *in vitro* and *in vivo*.<sup>10-12</sup>

### References

1. Brenner, S.L. and Korn, E.D. *J. Biol. Chem.* **255**(3), 841-844 (1980).
2. Lin, D.C., Tobin, K.D., Grumet, M., et al. *J. Cell Biol.* **84**, 455-460 (1980).
3. Ostlund, R.E., Jr., Leung, J.T., and Hajek, S.V. *J. Cell Biol.* **85**, 386-391 (1980).
4. Pinder, J.C. and Gratzner, W.B. *J. Cell Biol.* **96**(3), 768-775 (1983).
5. Flaumenhaft, R., Dilks, J.R., Rozenvayn, N., et al. *Blood* **105**(10), 3879-3887 (2005).
6. Taheri-Talesh, N., Horio, T., Araujo-Bazán, L., et al. *Mol. Biol. Cell* **19**, 1439-1449 (2008).
7. dos Santos, T., Varela, J., Lynch, I., et al. *PLoS One* **6**(9), 1-10 (2011).
8. Nightingale, T.D., White, I.J., Doyle, E.L., et al. *J. Cell Biol.* **194**(4), 613-629 (2011).
9. Walling, E.A., Krafft, G.A., and Ware, B.R. *Arch. Biochem. Biophys.* **264**(1), 321-332 (1988).
10. Yahara, I., Harada, F., Sekita, S., et al. *J. Cell Biol.* **92**, 69-78 (1982).
11. Chapla, V.M., Zeraik, M.L., Ximenes, V.F., et al. *Molecules* **19**, 6597-6608 (2014).
12. Deckert, M., Ticchioni, M., and Bernard, A. *J. Cell Biol.* **133**(4), 791-799 (1996).
13. Lee, J., Yi, J.-M., Kim, H., et al. *Biol. Pharm. Bull.* **37**(1), 6-12 (2014).

#### 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, 08/09/2023

#### 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