

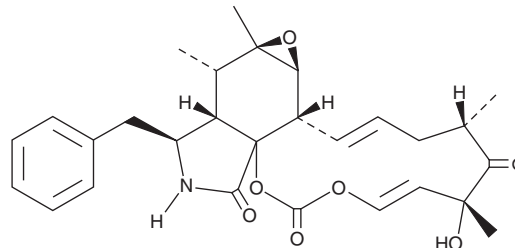
# PRODUCT INFORMATION



## Cytochalasin E

Item No. 11331

**CAS Registry No.:** 36011-19-5  
**Formal Name:** (1E,4S,6R,7E,11aS,14S,14aS,15S,15aR,16aS,16bS)-3,13,14,14a,15,15a,16a,16b-octahydro-6-hydroxy-4,6,15,15a-tetramethyl-14-(phenylmethyl)-[1,3]dioxacyclotridecino[4,5-d]oxireno[f]isoindole-5,10,12(4H,6H)-trione  
**Synonym:** NSC 175151  
**MF:** C<sub>28</sub>H<sub>33</sub>NO<sub>7</sub>  
**FW:** 495.6  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Aspergillus clavatus*



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

### Laboratory Procedures

Cytochalasin E is supplied as a crystalline solid. A stock solution may be made by dissolving the cytochalasin E in the solvent of choice, which should be purged with an inert gas. Cytochalasin E is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cytochalasin E in these solvents is approximately 2, 10, and 11 mg/ml, respectively.

Cytochalasin E is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cytochalasin E should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Cytochalasin E has a solubility of approximately 0.03 mg/ml in a 1:30 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

The cytochalasins are cell-permeable fungal metabolites which 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 E is an epoxide-containing analog of cytochalasin B (Item No. 11328) which potently and selectively inhibits the growth of endothelial cells (IC<sub>50</sub> < 1 nM), impairing angiogenesis and tumor growth.<sup>9</sup> This cytochalasin does not inhibit glucose transport or HIV-1 protease activity.<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**(2), 455-460 (1980).
3. Ostlund, R.E., Jr., Leung, J.T., and Hajek, S.V. *J. Cell Biol.* **85**(2), 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**(4), 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. Udagawa, T., Yuan, J., Panigrahy, D., et al. *J. Pharmacol. Exp. Ther.* **294**(2), 421-427 (2000).
10. Kletzien, R.F., Perdue, J.F., and Springer, A. *J. Biol. Chem.* **247**(9), 2964-2966 (1972).
11. Griffin, J.F., Rampal, A.L., and Jung, C.Y. *Proc. Natl. Acad. Sci. USA* **79**(12), 3759-3763 (1982).
12. Lombardo, Y.B., Drago, S., Chicco, A., et al. *Metabolism* **45**(12), 1527-1532 (1996).

#### 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, 08/11/2023

#### 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