

PRODUCT INFORMATION



Cytochalasin J

Item No. 18443

CAS Registry No.: 53760-20-6
Formal Name: 2,3,3a,4,5,6,6a,9,10,11,12,15-dodecahydro-6,12,15-trihydroxy-4,10,12-trimethyl-5-methylene-3-(phenylmethyl)-1H-cycloundec[d]isoindol-1-one

MF: C₂₈H₃₇NO₄

FW: 451.6

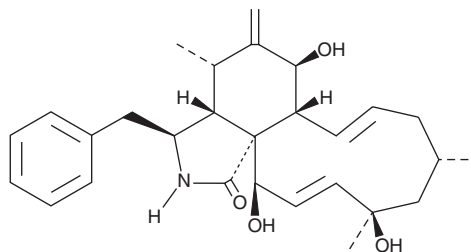
Purity: ≥95%

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 J is supplied as a lyophilisate. A stock solution may be made by dissolving the cytochalasin J in the solvent of choice, which should be purged with an inert gas. Cytochalasin J is soluble in organic solvents such as ethanol, methanol, DMSO and dimethyl formamide.

Description

The cytochalasins are cell-permeable fungal metabolites that inhibit actin polymerization.¹⁻⁴ This interferes with such diverse processes as cell movement, growth, phagocytosis, degranulation, and secretion.⁵⁻⁸ Cytochalasin J is a deacetylated analog of cytochalasin H that weakly inhibits actin assembly but significantly alters mitotic spindle microtubule organization and kinetochore structure.^{9,10} Like cytochalasin H, cytochalasin J suppresses the production of reactive oxygen species by stimulated human neutrophils.¹¹

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. *S 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. Wrench, G.A. and Snyder, J.A. *Cell Motil. Cytoskeleton* **36**(2), 112-124 (1997).
11. Chapla, V.M., Zeraik, M.L., Ximenes, V.F., et al. *Molecules* **19**, 6597-6608 (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