PRODUCT INFORMATION



Cytochalasin D

Item No. 11330

CAS Registry No.:	22144-77-0			
Formal Name:	(3S,3aR,4S,6S,6aR,7E,10S,12R,13E,15R,15			
	aR)-15-(acetyloxy)-3,3a,4,5,6,6a,9,10,12,15-			OH
	decahydro-6,12-dihydroxy-4,10,12-trimethyl-5-		[] [
	methylene-3-(phenylmethyl)-1H-cycloundec[d]	\sim		
	isoindole-1,11(2H)-dione		\checkmark \times	$\sim \sim \sim \uparrow^{\circ}$
Synonym:	NSC 209835			
MF:	C ₃₀ H ₃₇ NO ₆			
FW:	507.6		н о	OH
Purity:	≥95%			00
Supplied as:	A crystalline solid			Ť
Storage:	-20°C			
Stability:	≥2 years			
Information represents	s the product specifications. Batch specific analytical results	are provided or	n each certificate of	analysis

Laboratory Procedures

Cytochalasin D is supplied as a crystalline solid. A stock solution may be made by dissolving the cytochalasin D in the solvent of choice, which should be purged with an inert gas. Cytochalasin D is soluble in organic solvents such as DMSO and dichloromethane. The solubility of cytochalasin D in these solvents is approximately 100 and 10 mg/ml, respectively.

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 D is a cell-permeable inhibitor that binds actin filaments, but not actin monomers, to inhibit polymerization at concentrations as low as 0.2 μ M.² In this way, it prevents the migration of tumor cells.⁹

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 Gratzer, 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. Hayot, C., Debeir, O., Van Ham, P., et al. Toxicol. Appl. Pharmacol. 211, 30-40 (2006).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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.

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, 07/29/2021

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