


Cytochalasin D

Publication Number MAN0003709

Rev. 2.00








Catalog Number:	PHZ1063
Quantity:	1.0 mg
Lot Number:	See product label.
Appearance:	White solid
Molecular Formula:	C ₃₀ H ₃₇ NO ₆
Molecular Weight:	507.6
Purity:	98%
Summary:	Cytochalasin D is a cell permeable fungal toxin that binds to the barbed end of actin filaments inhibiting both the association and dissociation of subunits. This compound causes the disruption of actin filaments and inhibition of actin polymerization. Cytochalasin D is about 10–fold more effective than cytochalasin B, and does not inhibit monosaccharide transport across cell membranes.
Solubility:	Soluble in DMSO at a concentration of 25 mg/mL.
Sterility:	This product is not sterile.
Storage:	Store, as supplied, at –20°C. Upon solubilization, apportion into working aliquots and store at –20°C. Avoid repeated freeze/thaw cycles. Solutions are stable at –20°C for up to three months.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
References:	<p>Flanagan, M.D. and S. Lin (1980) Cytochalasins block actin filament elongation by binding to high affinity sites associated with F–actin. <i>J. Biol. Chem.</i> 255(3):835–838.</p> <p>Goddette, D.W. and C. Frieden (1986) Actin polymerization. The mechanism of action of cytochalasin D. <i>J. Biol. Chem.</i> 261(34):15974–15980.</p> <p>Cooper, J.A. (1987) Effects of cytochalasin and phalloidin on actin. <i>J. Cell Biol.</i> 105(4):1473–1478.</p> <p>Tangkijvanich, P., et al. (2002) Platelet-derived growth factor-BB and lysophosphatidic acid distinctly regulate hepatic myofibroblast migration through focal adhesion kinase. <i>Exp. Cell. Res.</i> 281(1):140–147.</p> <p>Goel, H.L., and C.S. Dey (2002) PKC-regulated myogenesis is associated with increased tyrosine phosphorylation of FAK, Cas, and paxillin, formation of Cas–CRK complex, and JNK activation. <i>Differentiation</i> 70(6):257–271.</p> <p>Tran, N.L., et al. (2002) Signal transduction from N–cadherin increases Bcl–2. Regulation of the phosphatidylinositol 3–kinase/Akt pathway by homophilic adhesion and actin cytoskeletal organization. <i>J. Biol. Chem.</i> 277(36):32905–32914.</p>
	Avoid contact with eyes, skin, and mucous membranes. Wear protective clothing when handling this product. Not for human use.







Related Products:

Product	Catalog no.
FAK Sample Pack	44-631
Akt/PKB [pS473] antibody	44-622G
JNK1&2 [pTpY183/185] antibody	44-682G
Paxillin [pY31] antibody	44-720G

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description
	Catalog Number
	Research Use Only
	Use by
	Manufacturer
	Without, does not contain
	Protect from light
	Directs the user to consult instructions for use (IFU), accompanying the product.

Symbol	Description
	Batch code
	In vitro diagnostic medical device
	Temperature limitation
	European Community authorized representative
	With, contains
	Consult accompanying documents

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

For Research Use Only. Caution: Not for human or animal therapeutic or diagnostic use.

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

Revision Date 13 May 2011

