

Cytochalasin D

Inhibitor of actin polymerization

Cytochalasin D is a cell-permeable mycotoxin that binds to the barbed end of actin filaments inhibiting both the association and dissociation of subunits. It causes the disruption of actin filaments and inhibition of actin polymerization. As a result, it impedes with a variety of processes including cell cycle, motility, phagocytosis, proliferation, amongst others. It is about 10-fold more effective than cytochalasin B (Prod. No. BML-T108). Cytochalasin D also inactivates low conductance K⁺ channels, modulates CD4 cross-linking in T lymphocytes, increases intracellular Ca²⁺ levels, and exhibits antitumor activity.

Citations: 52

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-T109-0001

1mg

Manuals, SDS & CofA

[View Online »](#)

- Highly pure and potent inhibitor of actin polymerization
- Blocker of low conductance potassium channel
- Highly cited

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NSC 209835, Zygosporin A
Appearance	White to off-white solid
CAS	22144-77-0
Couple Target	Potassium channel
Couple Type	Blocker
Formula	$C_{30}H_{37}NO_6$
MW	507.6
Purity	≥97% (HPLC, single spot by TLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol (5mg/ml), or methylene chloride (10mg/ml).
Source	Isolated from <i>Zygosporium mansonii</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com