

Cytochalasin A

HIV-1 protease inhibitor

Antibiotic. Fungal toxin. Inhibits glucose transport, actin polymerisation and blocks the formation of microtubuli. Inhibits cell division. Inhibits HIV-1 protease.

Citations: 5

[View Online »](#)

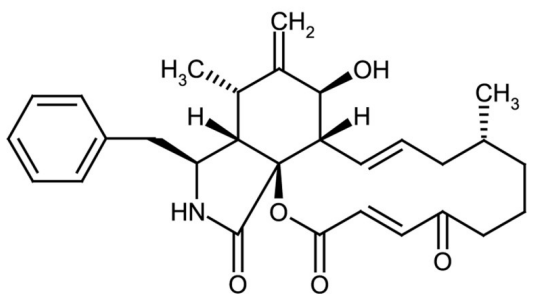
Ordering Information

[Order Online »](#)

ALX-380-057-M001	1mg
ALX-380-057-M010	10mg

Manuals, SDS & CofA

[View Online »](#)



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

Appearance	White to off-white powder.
CAS	14110-64-6
Couple Target	Protease
Couple Type	Inhibitor
Formula	$C_{29}H_{35}NO_5$
MI	14: 2790
MW	477.6
Purity	≥98% (HPLC, TLC)
Solubility	Soluble in acetone, DMSO or 100% ethanol.
Source	Isolated from <i>Drechslera dematoidea</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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