

Cytochalasin A

Mycotoxin produced by *Aspergillus*; actin polymerization inhibitor, Kv1.5 K⁺ channel blocker..

An anti-cytoskeletal drug which inhibits actin polymerization and has caused low stationary motility and membrane ruffling in K1735-M2 mouse melanoma cells.

Ordering Information

[Order Online »](#)

LKT-C9878-M001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dehydrophomin

CAS 14110-64-6

Couple Type Inhibitor

Formula $C_{29}H_{35}NO_5$

MW 477.6

Purity ≥98%



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