

Tentoxin

F-ATPase inhibitor

Cyclic tetrapeptide. Inhibitor of chloroplast F_1 -ATPase (CF_1). Does not inhibit the homologous enzymes from chloroplasts of insusceptible plant species or from bacteria or mitochondria.

Citations: 4

[View Online »](#)

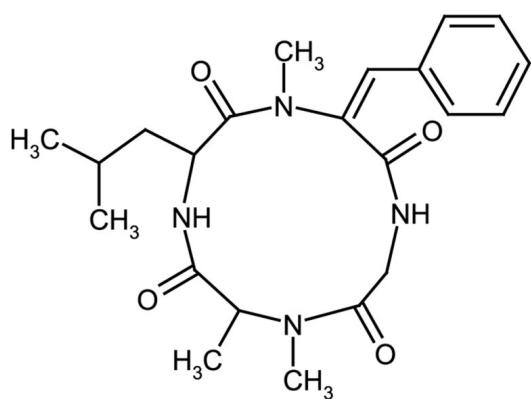
Ordering Information

[Order Online »](#)

ALX-350-132-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name cyclo-(N-Methyl-L-alanyl-L-leucyl-N-methyl-(Z)-dihydrophenylalanyl-glycyl)

Appearance Off-white to yellow powder.

CAS 28540-82-1

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{22}H_{30}N_4O_4$

MW 414.5

Purity ≥98% (HPLC, TLC)

Purity Detail Free of any peptide contaminations, especially dihydrotentoxin.

Solubility Soluble in DMSO (10mg/ml) or methanol (10mg/ml).

Source Isolated from *Alternaria alternata*.

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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