

Nonactin

Monovalent cation ionophore

Macrotetrolide antibiotic. Monovalent cation ionophore with high selectivity for ammonium and potassium. Specifically inhibits the processing of cytoplasmic precursor proteins destined for the mitochondria. Inhibits the effects on P170 glycoprotein-mediated efflux of chemotherapeutic agents in multidrug resistant (MDR) cancer cells.

Citations: 7

[View Online »](#)

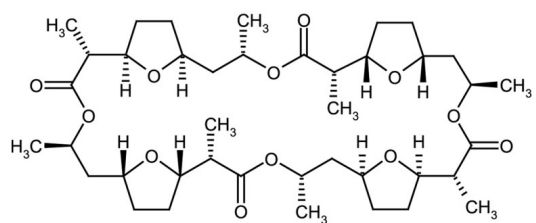
Ordering Information

[Order Online »](#)

ALX-450-008-M005	5mg
ALX-450-008-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Polynactin, Werramycin A
Appearance	White to off-white crystalline solid.
CAS	6833-84-7
Formula	$C_{40}H_{64}O_{12}$
MI	14: 6675
MW	736.9
Purity	≥98% (TLC) (Total homologs)
Purity Detail	Mixture of two homologs (~80% nonactin and ~20% monactin)
RTECS	RH1685000
Solubility	Soluble in DMSO (10mg/ml), dichloromethane (10mg/ml), methanol, 100% ethanol; insoluble in water.
Source	Isolated from <i>Streptomyces griseus</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-fr@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