

Prodigiosin

NADPH oxidase inhibitor

Antibacterial, antifungal, cytotoxic and antialgal activity. Has been shown to inhibit NADPH oxidase activation in macrophages. Induces apoptosis by phosphorylation of p38 MAP kinase. Additionally, inhibits the association of p47-phox and Rac (NADPH oxidase cytosolic components) to the plasma membrane in macrophages.

Citations: 6

[View Online »](#)

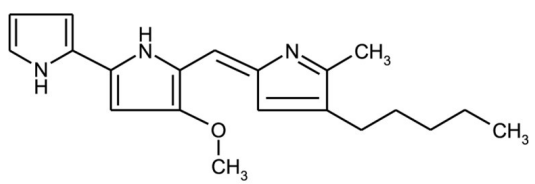
Ordering Information

[Order Online »](#)

ALX-380-321-MC25	0.25mg
ALX-380-321-M001	1mg

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

Alternative Name	Prodigiosine, 4-Methoxy-5-[(5-methyl-4-pentyl-2H-pyrrol-2-ylidene)methyl]-2,2'-bi[1H-pyrrole]
Appearance	Yellow to dark red solid.
CAS	82-89-3
Couple Target	NADPH oxidase
Couple Type	Inhibitor
Formula	$C_{20}H_{25}N_3O$
MI	14: 7767
MW	323.4
Purity	≥95% (HPLC)
Solubility	Soluble in dioxane, pyridine, methylene chloride, hexane, chloroform or methanol; insoluble in water.
Source	Isolated from <i>Serratia marcescens</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