

Molsidomine

NO releasing vasodilator

Long acting antianginal drug that is enzymatically converted in the liver to yield the active metabolite SIN-1 (Prod. No. BML-CN245). NO releasing vasodilator.

Citations: 6

[View Online »](#)

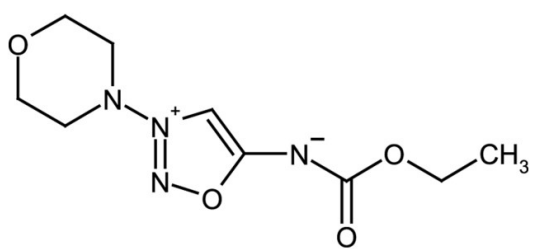
Ordering Information

[Order Online »](#)

ALX-430-003-M050	50mg
------------------	------

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	N-(Ethoxycarbonyl)-3-(4-morpholinyl)sydnoneimine, Morsidomine, SIN-10
Appearance	White solid.
CAS	25717-80-0
Formula	$C_9H_{14}N_4O_4$
Identity	Determined by NMR.
MI	14: 6234
MW	242.2
Purity	≥98% (TLC)
Solubility	Soluble in methanol.

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-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