

NOR-5

Nitric oxide donor

Nitric oxide (NO) donor. The half-life time of 0.5mM NOR-5 is 20 hours at 37°C in 0.1M PBS, pH 7.4.

Citations: 4

[View Online »](#)

Ordering Information

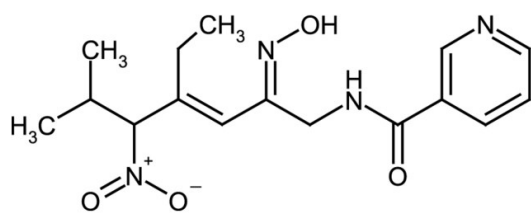
[Order Online »](#)

ALX-430-125-M010

10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-2-((E)-4-Ethyl-3[(Z)-hydroxyimino]6-methyl-5-nitro-heptenyl)-3-pyridinecarboxamide
Appearance	White to off-white powder.
Formula	$C_{16}H_{22}N_4O_4$
MW	334.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).

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