

# NOR-3

## Nitric oxide donor

Non-thiol-based nitric oxide (NO) donor. NOR-3 shows a variety of biological and pharmacological effects. Strong vasodilatory effects in rat aorta, rabbit aorta and dog coronary artery; potent inhibitory activities against antiplatelet aggregation and thrombus formation; effective aggregation inhibition of human platelet raising by ADP; cardioprotective effects in the ischemia/reperfusion system; antianginal effect.

Citations: 16

[View Online »](#)

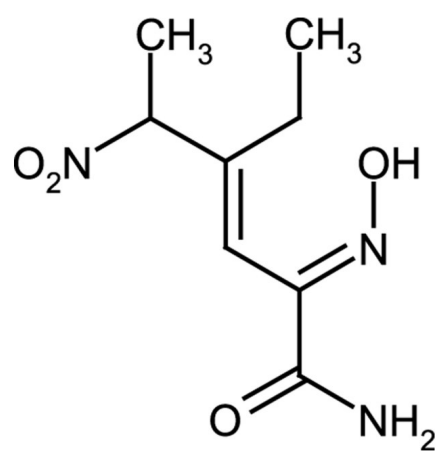
## Ordering Information

[Order Online »](#)

ALX-430-011-M005	5mg
ALX-430-011-M010	10mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Crystals are sensitive to moisture and become discolored on exposure to air. Solutions are stable for 1 month when stored at -20°C. Solvents should be purged with inert gas prior to addition of compound.
<b>Handling</b>	Protect from light and moisture. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	FK409, (±)-(E)-4-Ethyl-2-[(E)-hydroxyimino]-5-nitro-3-hexeneamide
<b>Appearance</b>	White to off-white crystalline solid.
<b>CAS</b>	138472-01-2
<b>Formula</b>	$C_8H_{13}N_3O_4$
<b>Identity</b>	Identity determined by NMR.
<b>MW</b>	215.2
<b>Purity</b>	≥97% (HPLC)
<b>Solubility</b>	Soluble in DMSO (10mg/ml).



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