

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 »](#)

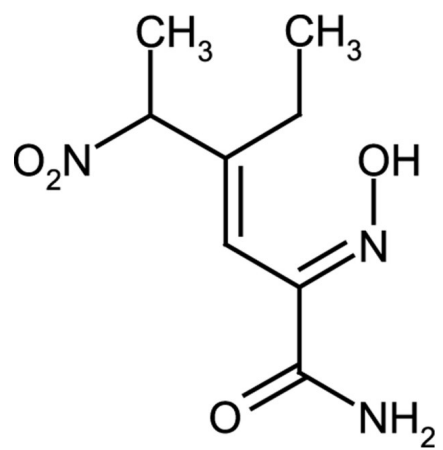
Ordering Information

[Order Online »](#)

ALX-430-011-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Handling	Protect from light and moisture. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	FK409, (±)-(E)-4-Ethyl-2-[(E)-hydroxyimino]-5-nitro-3-hexeneamide
Appearance	White to off-white crystalline solid.
CAS	138472-01-2
Formula	$C_8H_{13}N_3O_4$
Identity	Identity determined by NMR.
MW	215.2
Purity	≥97% (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-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