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 »

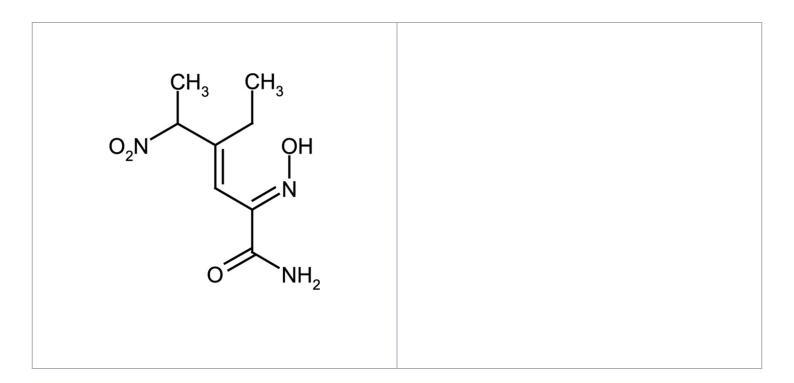
Ordering Information

Order Online »

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

Manuals, SDS & CofA

View Online »



Handling & Storage

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)

Soluble in DMSO (10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 infouk@enzolifesciences.com