

DEA NONOate

Nitric oxide donor

Nitric oxide (NO) donor.

Citations: 22

[View Online »](#)

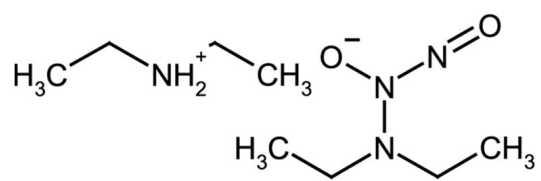
Ordering Information

[Order Online »](#)

ALX-430-034-M010	10mg
ALX-430-034-5005	5x5mg
ALX-430-034-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year. Aqueous solutions should be prepared fresh daily.
Handling	Protect from light and oxygen. Packaged under inert gas. Hygroscopic.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DEA/NO, 2-(N,N-Diethylamino)-diazene-2-oxide . diethylammonium salt
Appearance	White to off-white solid.
CAS	372965-00-9
Formula	$C_4H_{10}N_3O_2 \cdot C_4H_{12}N$
MW	132.1 . 74.1
Purity	≥97% (NMR)
Solubility	Soluble in 100% ethanol (25mg/ml), DMSO (2mg/ml), dimethyl formamide (2mg/ml) or PBS, pH 7.2 (10mg/ml).

Technical Info / Product Notes

Dissociates to the free amine and nitric oxide in a pH-dependent manner following first order kinetics. To initiate the release of nitric oxide, add stock alkaline solution of DEA NONOate to excess buffer of pH 7.0 – 7.4. The half life in 0.1 M phosphate buffer, pH 7.4, is $t_{1/2} = 16$ min. at 22 °C or $t_{1/2} = 2$ min. at 37 °C. Liberates 1.5 mol of NO per mol of parent compound. Decomposition is nearly instantaneous at pH 5.0.



Last modified: May 29, 2024
usa@enzolifesciences.com

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@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