

PAPA NONOate

Nitric oxide donor

Nitric oxide (NO) donor. The half-life time of PAPA NONOate is 76.6 min. at 22°C in 0.1M phosphate buffer, pH 7.4. Decomposition of NONOates is nearly instantaneous at pH 5.0.

Citations: 8

[View Online »](#)

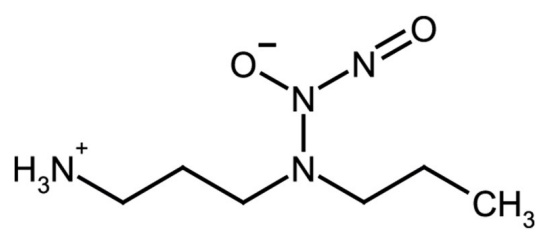
Ordering Information

[Order Online »](#)

ALX-430-016-M005	5mg
ALX-430-016-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Relatively stable in alkaline solution.
Handling	Protect from moisture. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PAPA/NO, NOC-15, (Z)-1-[N-(3-Ammoniopropyl)-N-(n-propyl)amino]diazene-1-ium-1,2-diolate
Appearance	White to off-white solid.
CAS	146672-58-4
Formula	$C_6H_{16}N_4O_2$
MW	176.2
Purity	≥97% (TLC)
Solubility	Soluble in water or methanol.

Last modified: May 29, 2024



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