

DPTA NONOate

Nitric oxide donor

Nitric oxide (NO) donor. The half-life time is 5 hours at 22°C in 0.1M phosphate buffer, pH 7.4. Decomposition of NONOates is nearly instantaneous at pH 5.0.

Citations: 5

[View Online »](#)

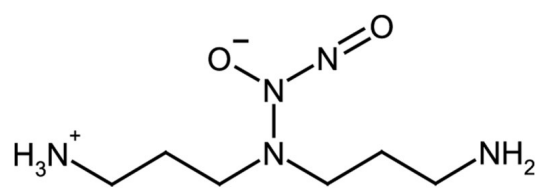
Ordering Information

[Order Online »](#)

ALX-430-066-M005	5mg
ALX-430-066-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture. Aqueous solutions should be prepared fresh daily.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dipropylenetriamine NONOate
Appearance	White to off-white solid.
CAS	146724-95-0
Formula	$C_6H_{17}N_5O_2$
MW	191.2
Purity	≥98% (NMR)
Solubility	Highly soluble in water.

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)