

# DETA NONOate

## Nitric oxide donor

Nitric oxide (NO) donor. Induces apoptosis in macrophages. The half-life time of DETA NONOate is 20 hours at 37°C and 56 hours at 22-25°C in 0.1M phosphate buffer, pH 7.4. Decomposition of NONOates is nearly instantaneous at pH 5.0. DETA NONOate liberates 2 mol of NO per mol of parent compound.

Citations: 24

[View Online »](#)

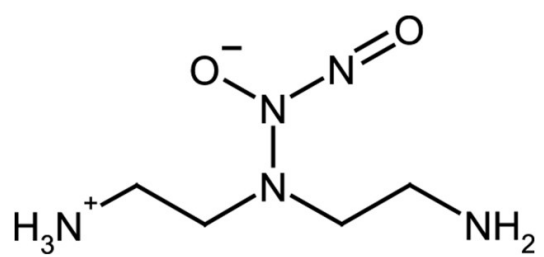
## Ordering Information

[Order Online »](#)

|                  |       |
|------------------|-------|
| ALX-430-014-M005 | 5mg   |
| ALX-430-014-5005 | 5x5mg |
| ALX-430-014-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.                                 |
| Long Term Storage | -80°C                                                            |
| Shipping          | Dry Ice                                                          |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                                                                                                                                          |
|------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Alternative Name | NOC-18, DETA/NO, (Z)-1-[2-(2-Aminoethyl)-N-(2-ammonioethyl)amino]diazene-1-ium-1,2-diolate, 3,3-Bis(aminoethyl)-1-hydroxy-2-oxo-1-triazene, 2,2'-(Hydroxynitrosohydrazino)bis-ethanamine |
| Appearance       | White to off-white solid.                                                                                                                                                                |
| CAS              | 146724-94-9                                                                                                                                                                              |
| Formula          | $C_4H_{13}N_5O_2$                                                                                                                                                                        |
| MW               | 163.2                                                                                                                                                                                    |
| Purity           | ≥97% (1H-NMR)                                                                                                                                                                            |
| RTECS            | KH2502500                                                                                                                                                                                |
| Solubility       | Soluble in water or ethanol.                                                                                                                                                             |

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