

Spermine NONOate

Nitric oxide donor

Nitric oxide (NO) donor. A convenient reagent for preparing aqueous solutions of NO.

Citations: 13

[View Online »](#)

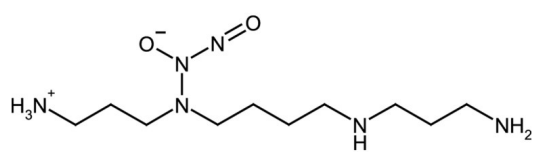
Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-430-013-M005 | 5mg |
| ALX-430-013-5005 | 5x5mg |
| ALX-430-013-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 light and oxygen. Keep under inert gas. Hygroscopic. |
| Long Term Storage | -80°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|--|
| Alternative Name | SPER/NO, (Z)-1-{N-[3-Aminopropyl]-N-[4-(3-aminopropylammonio)butyl]-amino}-diazene-1-ium-1,2-diolate |
| Appearance | White to off-white crystalline solid. |
| CAS | 136587-13-8 |
| Formula | $C_{10}H_{26}N_6O_2$ |
| MW | 262.4 |
| Purity | ≥97% (NMR) |
| Solubility | Soluble in water, methanol or 100% ethanol. |

Technical Info / Product Notes

Product dissociates to the free amine and nitric oxide in a pH dependent manner following first order kinetics. Alkaline solutions of NONOates (in 0.01M NaOH) are stable and can be stored at 0°C for 24 hours. To initiate the release of nitric oxide add alkaline solution of Spermine NONOate to excess buffer of pH 7.0 – 7.4. Relatively high concentrated solutions can be prepared for further dilution. Half life in 0.1M phosphate buffer (pH 7.4) at 37 C is $t_{1/2}$ (min.) 39. Decomposition is nearly instantaneous at pH 5.



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

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