Spermine NONOate

Nitric oxide donor

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

Citations: 13

View Online »

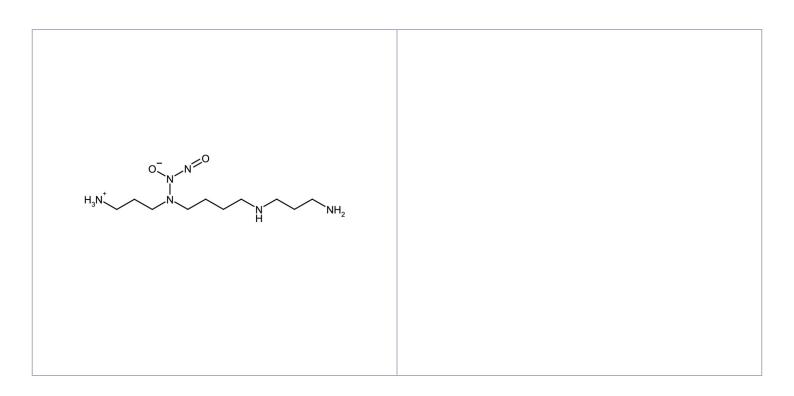
Ordering Information

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}-diazen-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 NotesProduct 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 t½ (min.) 39. Decomposition is nearly

instantaneous at pH 5.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com