

NOC-7

Nitric oxide donor

Nitric oxide (NO) donor. The half-life time of 0.1mM NOC-7 is 5 min. at 37°C in 0.1M PBS, pH 7.4.

Citations: 3

[View Online »](#)

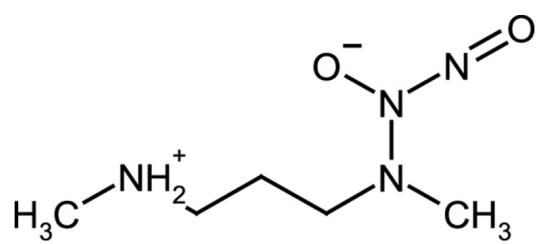
Ordering Information

[Order Online »](#)

ALX-430-018-M005	5mg
ALX-430-018-46190	5mg
ALX-430-018-M010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare fresh stock solution prior to use. Store in ice bath during the experiment. Stock solution will lose 5% of activity per day when stored at -20°C.
Handling	Protect from light and moisture. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	1-Hydroxy-2-oxo-3-(N-3-methyl-aminopropyl)-3-methyl-1-triazene
Appearance	White to off-white crystalline solid.
CAS	146724-84-7
Formula	$C_5H_{14}N_4O_2$
MW	162.2
Purity	≥96% (HPLC)
Reconstitution	Reconstitute with 0.1M sodium hydroxide solution to 10-50mM.

Last modified: May 29, 2024



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