

Diethyldithiocarbamic acid . sodium salt . trihydrate (high purity)

NO spin trapping reagent

Nitric oxide (NO) spin-trapping reagent. Thiol and iron chelator. Inhibits induction of macrophage nitric oxide synthase (NOS). Has been shown to be an inhibitor of the nuclear transcription factor κ B (NF- κ B).

Citations: 17

[View Online »](#)

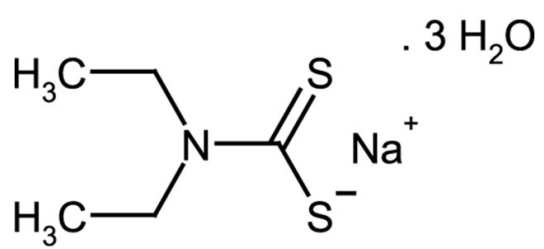
Ordering Information

[Order Online »](#)

ALX-400-003-G025	25g
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DETC . 3H₂O (high purity), Diethyldithiocarbamate . 3H₂O (high purity)

Appearance White to off-white solid.

CAS 20624-25-3

Couple Target NF-kappaB

Couple Type Inhibitor

Formula C₅H₁₀NS₂ . Na . 3H₂O

MI 14: 3378

MW 148.3 . 23.0 . 54.0

Purity ≥97% (Titration)

RTECS EZ6550000

Solubility Soluble in water (20mg/ml).

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-be@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)