

MGD . sodium salt . monohydrate

Spin trapping reagent

Together with FeSO_4 MGD is a useful component for the formation of the $\text{MGD}_2\text{-Fe}^{2+}$ complex, which is an excellent nitric oxide (NO) spin-trapping reagent. The $\text{MGD}_2\text{-Fe}^{2+}$ complex is quite unstable, especially in the presence of dissolved oxygen. Thus, the complex should be used immediately after being made. An excess (usually 5-fold excess), of MGD to Fe^{2+} is used for making the complex with FeSO_4 to give a more stable complex solution. Acidic conditions should be avoided because dithiocarbamate tends to decompose forming toxic carbon disulfide. It was reported that MGD and $\text{Fe}(\text{MGD})_2$ do not exhibit toxicity up to 8mmol/kg and 0.3mmol/kg, respectively.

Citations: 15

[View Online »](#)

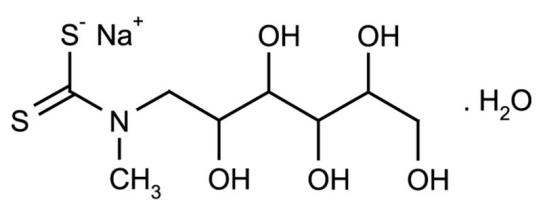
Ordering Information

[Order Online »](#)

ALX-400-014-M050	50mg
ALX-400-014-M250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(Dithiocarbamoyl)-N-methyl-D-glucamine . Na . H ₂ O, N-Methyl-D-glucamine dithiocarbamate
Appearance	White to off-white solid.
CAS	91840-27-6
Formula	C ₈ H ₁₆ NO ₅ S ₂ Na . H ₂ O
MW	293.3 . 18.0
Purity	≥98% (HPLC)
Solubility	Soluble in water.

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)