

GSNO

Nitric oxide donor

NO donor which spontaneously releases NO under physiological conditions. Half-life of 159 hours under physiological conditions. Selectively inhibits platelet activation and displays cardioprotective effects. Inhibits NF- κ B activation. Inhibits endothelial cell proliferation and induces apoptosis in T cells.

Citations: 25

[View Online »](#)

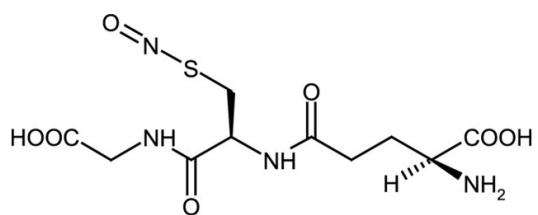
Ordering Information

[Order Online »](#)

ALX-420-002-M025	25mg
ALX-420-002-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name S-Nitroso-L-glutathione, N-(N-L-γ-Glutamyl-S-nitroso-L-cysteinyl)-glycine

Appearance Crystalline solid.

CAS 57564-91-7

Formula $C_{10}H_{16}N_4O_7S$

MW 336.3

Purity ≥98% (HPLC)

Solubility Soluble in water (25mg/ml), DMSO (20mg/ml) or PBS; sparingly soluble in 100% ethanol, methanol or dimethyl formamide.

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