

RSSR

Supersensitive spin label

Cell permeable supersensitive spin label (detection limit varies from 10nM to 100nM) for quick detection of reduced thiols using ESR. Allows following the reactions of sulfhydryl groups with RSSR to form thiol spin label adducts, for the monitoring of intracellular redox states of glutathione and other thiols.

Citations: 13

[View Online »](#)

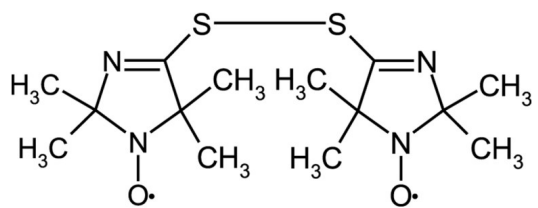
Ordering Information

[Order Online »](#)

ALX-430-102-M025	25mg
------------------	------

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	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	bis-(2,2,5,5-Tetramethyl-3-imidazoline-1-oxyl-4-yl)disulfide, biradical
Appearance	Off-white to yellow crystalline powder.
CAS	123557-49-3
Formula	$C_{14}H_{24}N_4O_2S_2$
MW	344.5
Purity	≥97% (HPLC, GLC)
Solubility	Soluble in DMSO, water, 100% ethanol or methanol; DMSO is recommended for stock solution.



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-
be@enzolifesciences.com](mailto:info-be@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)