

MTSSL

Thiol-specific spin label

Highly reactive thiol-specific spin label. Has been used to label cysteine residues in proteins (site-directed labeling, SDS-labeling). Allows protein structure and protein dynamics determination as well as the study of protein-protein and protein-oligonucleotide interactions.

Citations: 24

[View Online »](#)

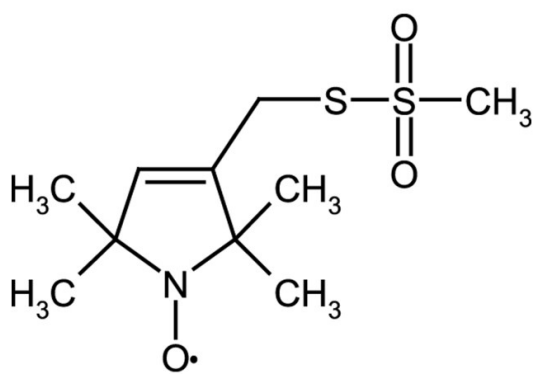
Ordering Information

[Order Online »](#)

ALX-430-134-M010	10mg
ALX-430-134-M050	50mg

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(1-Oxyl-2,2,5,5-tetramethylpyrroline-3-methyl) methanethiosulfonate
Appearance	Yellow crystalline solid.
CAS	81213-52-7
Formula	$C_{10}H_{18}NO_3S_2$
MW	264.3
Purity	≥98% (HPLC)
Solubility	Soluble in water, methanol, 100% ethanol, DMSO, acetonitrile or acetone.

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-be@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