

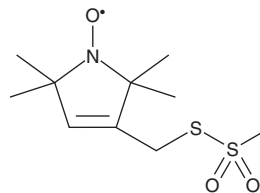
PRODUCT INFORMATION



MTSSL

Item No. 16463

CAS Registry No.: 81213-52-7
Formal Name: 2,5-dihydro-2,2,5,5-tetramethyl-3-
[[[(methylsulfonyl)thio]methyl]-1H-pyrrol-1-yloxy
MF: C₁₀H₁₈NO₃S₂
FW: 264.4
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MTSSL is supplied as a crystalline solid. A stock solution may be made by dissolving the MTSSL in the solvent of choice, which should be purged with an inert gas. MTSSL is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MTSSL in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

MTSSL is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MTSSL should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MTSSL has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MTSSL is a paramagnetic nitroxide spin label featuring a methanethiosulfonate group for site-directed spin labeling, particularly through disulfide linkages with cysteine residues on proteins.^{1,2} When combined with electron paramagnetic resonance spectroscopy, site-directed spin labeling is used to study the structural dynamics of proteins, including membrane proteins.^{1,3}

References

1. Perozo, E., Cortes, D.M., and Cuello, L.G. Structural rearrangements underlying K⁺-channel activation gating. *Science* **285**(5424), 73-78 (1999).
2. Gaponenko, V., Howarth, J.W., Columbus, L., *et al.* Protein global fold determination using site-directed spin and isotope labeling. *Protein Sci.* **9**(2), 302-309 (2000).
3. Tombolato, F., Ferrarini, A., and Freed, J.H. Dynamics of the nitroxide side chain in spin-labeled proteins. *J. Phys. Chem. B* **110**(51), 26248-26259 (2006).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/09/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM