

ReadiUse™ Traut's Reagent (2-iminothiolane) *100 mM DMSO Solution*

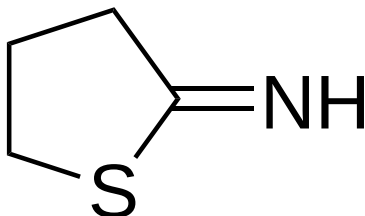
Catalog Number: 4514

Unit Size: 5 mL

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	137.63
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

ReadiUse™ Traut's Reagent (2-iminothiolane) is a cyclic thioimide thiolation reagent supplied as a 100 mM solution in DMSO for introducing sulfhydryl groups onto biomolecules containing primary amines. The reagent reacts selectively with lysine side chains and N-terminal amino groups in aqueous buffers near neutral pH (pH7-10) to generate thiol-functionalized biomolecules. During modification, the reagent preserves charge properties similar to the original amino group, helping maintain biomolecule solubility and functionality.

The introduced sulfhydryl groups can subsequently participate in thiol-reactive conjugation chemistries including maleimide-based crosslinking reactions. ReadiUse™ Traut's Reagent can be used for thiolation of antibodies, proteins, peptides, and other amino-containing biomolecules in bioconjugation, labeling, immobilization, and crosslinking workflows. The reagent introduces a short spacer arm that can help reduce steric hindrance during downstream conjugation reactions.