

## SATA (N-succinimidyl S-acetylthioacetate)

Catalog Number: 4515

Unit Size: 25 mg

### Product Details

---

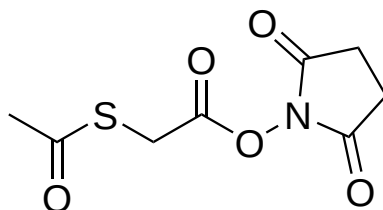
Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	231.23
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

ReadiUse™ SATA (N-succinimidyl S-acetylthioacetate) is an NHS ester-based thiolation reagent used for introducing protected sulfhydryl groups onto proteins and other primary amine-containing biomolecules. The NHS ester reacts with lysine side chains and N-terminal amino groups to form stable covalent amide bonds while introducing an acetyl-protected sulfhydryl functionality. The protected thiol group can subsequently be deprotected using hydroxylamine to generate a free sulfhydryl group for downstream conjugation reactions.

The protected sulfhydryl strategy enables preparation and storage of modified biomolecules prior to thiol exposure and conjugation. ReadiUse™ SATA can be used for thiolation of antibodies, proteins, peptides, and other amino-containing biomolecules in crosslinking, labeling, immobilization, and bioconjugation workflows. Following deprotection, the exposed sulfhydryl groups can participate in thiol-reactive chemistries including maleimide-based conjugation and disulfide bond formation.