

## Acetic Acid N-Succinimidyl Ester

Catalog Number: 4508

Unit Size: 100 mg

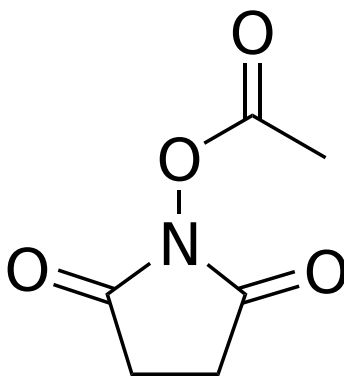
### Product Details

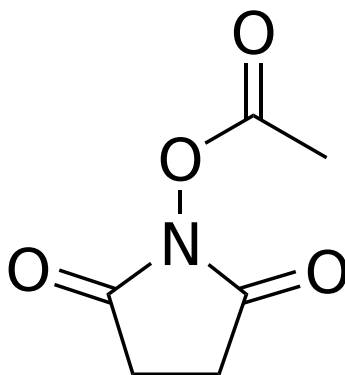
---

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

### Chemical Properties

---

Appearance	Solid
Molecular Weight	157.13
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

Acetic Acid N-Succinimidyl Ester is a convenient reagent for blocking amino groups of biological molecules. It has an extremely small size, thus minimizing the negative effect of the molecules to be modified. It readily reacts with biological molecules (such as antibodies, peptides, amino-modified oligos) under mild conditions with optimized pH from 7.5 to 9.0.