

## 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-KLH Conjugate

Catalog number: 50014

Unit size: 1 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Liquid
------------	--------

### Applications

---

3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-KLH Conjugate is used for developing anti-3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ) antibodies that are typically used in an ELISA assay. 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-BSA Conjugate has multiple 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ) molecules on each KLH to maximize its immunogenicity.