

3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-HRP Conjugate

Catalog Number: 50015

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	N/A

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-HRP conjugate is used for developing an ELISA assay for quantifying 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ). 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ)-HRP conjugate has minimal number of 3-Amino-5-morpholinomethyl-2-oxazolidinone (AMOZ) molecules on each HRP to maximize its assay sensitivity.