

3-Amino-2-oxazolidine (AOZ)-BSA Conjugate

Catalog number: 50004
Unit size: 1 mg

Product Details

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

Chemical Properties

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

Applications

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