

Cinoxacin-BSA Conjugate

Catalog number: 50023

Unit size: 1 mg

Product Details

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Liquid

Applications

Cinoxacin-BSA Conjugate is used for developing anti-cinoxacin antibodies that are typically used in an ELISA assay. Cinoxacin-BSA Conjugate has multiple cinoxacin molecules on each BSA to maximize its immunogenicity.