

Cinoxacin-KLH Conjugate

Catalog number: 50024
Unit size: 1 mg

Product Details

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

Chemical Properties

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

Applications

Cinoxacin-KLH 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 KLH to maximize its immunogenicity.