

## **Spectinomycin-KLH Conjugate**

Catalog number: 50039

Unit size: 1 mg

**Product Details** 

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Liquid

## **Applications**

Spectinomycin-KLH conjugate is used for developing anti-spectinomycin antibodies that are typically used in an ELISA assay. Spectinomycin-KLH conjugate has multiple spectinomycin molecules on each BSA to maximize its immunogenicity.