

Sulfathiazol-KLH Conjugate

Catalog number: 50030

Unit size: 1 mg

Product Details

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Liquid

Applications

Sulfathiazol-KLH conjugate is used for developing anti-sulfathiazol antibodies that are typically used in an ELISA assay. The anti-sulfathiazol antibodies may also recognize sulfatroxazol, sulfamethoxazol, sulfamethiazol and sulfisoxazole. Sulfathiazol-KLH conjugate has multiple sulfathiazol molecules on each KLH to maximize its immunogenicity.