

Florfenicol-BSA Conjugate

Catalog number: 50035

Unit size: 1 mg

Product Details

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Liquid

Applications

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