

Chloramphenicol-Ovalbumin Conjugate

Catalog number: 50000

Unit size: 1 mg

Product Details

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

Chemical Properties

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

Applications

Chloramphenicol-Ovalbumin Conjugate is used for developing anti-chloramphenicol antibodies that are typically used in an ELISA assay. Chloramphenicol-Ovalbumin Conjugate has multiple chloramphenicol molecules on each ovalbumin to maximize its immunogenicity.