

Anti-Enrofloxacin antibody *Mouse anti-human, monoclonal IgG2b*

Catalog number: V100260, V100261

Unit size: 50 ug, 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Soluble In	Water

Applications

This antibody was prepared in mouse with reactivity against human. Applications include ELISA.