

Anti-Clenbuterol antibody *Mouse anti-human, monoclonal IgG1*Catalog number: V100245, V100246
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.