

**Anti-TIMP1 antibody *Mouse anti-human,
monoclonal IgG2a***Catalog number: V100070, V100071
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, CLIA.