

**Anti-Furazolidone antibody \*Mouse anti-human, monoclonal IgG1\***Catalog number: V100250, V100251  
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.