

**Purified Mouse Anti-dog Intra-Acrosomal Protein Antibody \*Ds-2, monoclonal\***

Catalog number: V1031575

Unit size: 0.1 mg

**Product Details**

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Dog
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Intra-Acrosomal Protein
Clone	Ds-2

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ICC, WB

**Applications**

---

Acrosomal protein SP-10 is a protein with a molecular weight of 33 kDa, found in the acrosomal vesicle of cells. In humans, acrosomal protein SP-10 plays an important role in organismal processes, in particular, multicellular organism development.