

**Purified Mouse Anti-human NHERF1  
Antibody \*EBP-10, monoclonal\***

Catalog number: V1031850

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	NHERF1
Clone	EBP-10

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IP, WB, IHC(P)

**Applications**

---

This is an antibody that targets NHERF1.