

**Purified Mouse Anti-mammalian  
Neurofilament H Antibody \*NF-01,  
monoclonal\***Catalog number: V1031825  
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	Mammalian
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Neurofilament H
Clone	NF-01

**Biological Properties**

---

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

**Applications**

---

This is an antibody that targets Neurofilament H.