

**Purified Mouse Anti-human/rat/pig
Neurofilament H Antibody *NF-05,
monoclonal, Cross Adsorbed***

Catalog number: V1031830
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, rat, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Neurofilament H
Clone	NF-05

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with under optimal conditions
Application	WB, IHC(F), ELISA

Applications

This is an antibody that targets Neurofilament H.