

**Purified Mouse Anti-human/mouse/pig
SLP76 Antibody *SLP-76/03, monoclonal***

Catalog number: V1032125

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, mouse, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	SLP76
Clone	SLP-76/03

Biological Properties

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

Applications

This is an antibody that targets SLP76.