

Purified Rabbit Anti-mouse Blnk Antibody
PAb (411), monoclonalCatalog number: V103255
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rabbit
Immunogen	Blnk
Clone	PAb (411)

Biological Properties

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

Applications

B-cell linker protein is a protein expressed in the cytoplasmic ribonucleoprotein granule, plasma membrane and cytoplasm of cells. Sequencing of B-cell linker protein has shown it contains a primary structural unit, the SH2 domain. It has been found to be involved in organismal processes, for example, intracellular signal transduction and B cell activation. B-cell linker protein is the subject of comprehensive application because of the fact that it is a component of the transmembrane receptor protein tyrosine kinase signaling pathway.