

**Purified Mouse Anti-species independent
Phosphotyrosine Antibody *P-Tyr-01,
monoclonal***Catalog number: V1032010
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	Species independent
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Phosphotyrosine
Clone	P-Tyr-01

Biological Properties

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

Applications

O-Phospho-L-tyrosine (phospho-L-tyrosine, phosphotyrosine) is a phosphorylated amino acid that occurs in a number of proteins. Phospho-L-tyrosine is vital to the research of important biological receptors like cell-surface receptor, particularly in cell-lines such as hematopoietic cells. Furthermore, it is vital to the application of vital cellular processes like protein synthesis and it is essential to the application of key cell-based activities like SRC kinase activity and kinase activity.