

Purified Mouse Anti-human Vimentin Antibody *VI-RE/1, monoclonal, Cross Adsorbed*

Catalog number: V1032435

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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Vimentin
Clone	VI-RE/1

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse, pig serum and then conjugated with under optimal conditions
Application	FC (QC TESTED), WB, ICC, ELISA

Applications

Vimentin is a 54 kDa protein that can be found in the cytosol, extracellular exosome and cytoskeleton of cells. In Homo sapiens, vimentin binds to double-stranded RNA and identical protein. It is the subject of intensive research in part because it is a member of the cytokine-mediated signaling pathway. Vimentin is thought to be essential to organismal processes, for example, cellular response to interferon- γ , SMAD protein signal transduction and astrocyte development. Sequencing of vimentin has supported it contains a primary structural unit, the IF rod domain. Vimentin is thought to be essential to mRNA stability. It is an inhibitor of neuron projection development but also is a promoter of translation and collagen biosynthetic process. Vimentin is clinically significant because abnormalities in its function have been thought to be involved with diseases like Cataract 30, multiple types (CTRCT30).