

**Purified Mouse Anti-human/mouse/pig
Vimentin Antibody *VI-10, monoclonal***Catalog number: V1032440
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse, pig, chicken, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Vimentin
Clone	VI-10

Biological Properties

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

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).