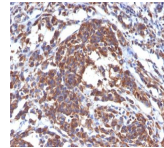


# Vimentin Monoclonal Mouse Antibody (LN-6)



## Product Description

This MAb reacts with a 58 kDa protein identified as vimentin. It reacts with a non-hematopoietic epitope of vimentin and shows no cross-reaction with other closely related intermediate filament proteins (IFP's) such as desmin, keratin, neurofilament, and glial fibrillary acid protein. Vimentin is ubiquitously expressed in mesenchymal cells such as fibroblasts, smooth muscle cells, and endothelium. Antibody against vimentin is useful as part of an antibody panel for differential diagnosis of tumors of unknown origin. Ab-2 does not react with leukocyte common antigen-positive tissues such as lymphomas and leukemias.

This antibody is available purified, with BSA (0.2 mg/mL) or purified, BSA-free (1 mg/mL).

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Histochem Cytochem (1989 ) 37(9) 1363-1370. (IHC, FFPE; IHC, frozen; IF; IP)
2. Int J Toxicol (2007) 26:289-295. (IHC; WB)
3. Stem Cell Res Ther (2018) 9: 178 (Flow, intracellular)

## Product attributes

<b>Antibody number</b>	#0060
<b>Antibody reactivity (target)</b>	Vimentin
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	LN-6
<b>Isotype</b>	IgM
<b>Molecular weight</b>	57-60 kDa
<b>Synonyms</b>	VIM
<b>Human gene symbol</b>	VIM
<b>Entrez gene ID</b>	7431
<b>SwissProt</b>	P08670
<b>Unigene</b>	455493
<b>Immunogen</b>	Human thymic nuclear extract
<b>Antibody target cellular fractionation</b>	Cytoskeleton
<b>Species reactivity</b>	Cat, Cow, Human, Mouse, Pig, Rabbit, Rat, Sheep
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Expected antibody applications</b>	IHC (frozen) (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
<b>Antibody application notes</b>	Immunohistology formalin-fixed 1-2 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	Jurkat cells, Sarcomas, Melanomas
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody research areas</b>	Cancer, Cytoskeleton
<b>Antibody/conjugate formulation</b>	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide