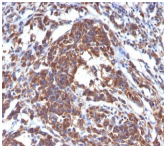


Vimentin Monoclonal Mouse Antibody (LN-6)



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

Antibody number	#0060
Antibody reactivity (target)	Vimentin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LN-6
Isotype	IgM
Molecular weight	57-60 kDa
Synonyms	VIM
Human gene symbol	VIM
Entrez gene ID	7431
SwissProt	P08670
Unigene	455493
Immunogen	Human thymic nuclear extract
Antibody target cellular localization	Cytoskeleton
Species reactivity	Cat, Cow, Human, Mouse, Pig, Rabbit, Rat, Sheep
Verified antibody applications	IHC (FFPE) (verified)
Expected antibody applications	IHC (frozen) (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
Antibody application notes	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
Positive control	Jurkat cells, Sarcomas, Melanomas
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody research areas	Cancer, Cytoskeleton
Antibody/conjugate formulation	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