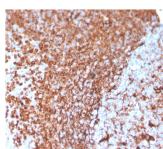


Vimentin Monoclonal Mouse Antibody (VIM/3736)



Product Description

This MAbs reacts with a 58 kDa protein identified as vimentin. It shows no cross-reaction with other closely related intermediate filament proteins (IFP s) such as desmin, keratin, neurofilament, and glial fibrillary acid protein. Anti-vimentin alone is of limited value as a diagnostic tool; however, when used in panels with other antibodies, it is useful for the sub-classification of a given tumor. Expression of vimentin, when used in conjunction with anti-keratin, is helpful when distinguishing melanomas from undifferentiated carcinomas and large cell lymphomas. All melanomas and Schwannomas react strongly with anti-vimentin. It labels a variety of mesenchymal cells, including melanocytes, lymphocytes, endothelial cells, and fibroblasts. Non-reactivity of anti-vimentin is often considered more useful than its positive reactivity, since there are a few tumors that do not contain vimentin, e. g. hepatoma and seminoma. Anti-vimentin is also useful as a tissue process control reagent.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3736, Anti-Vimentin (VIM/3736)

Product attributes

Antibody number	#3736
Antibody reactivity (target)	Vimentin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	VIM/3736
Isotype	IgG2b, kappa
Molecular weight	57-60 kDa
Synonyms	VIM
Human gene symbol	VIM
Entrez gene ID	7431
SwissProt	P08670
Unigene	455493
Immunogen	A recombinant fragment (around aa 2-466) of human Vimentin protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human
Positive control	U87, A375, Raji, HEP2 cells. Sarcomas or Melanomas.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
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
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on September 2, 2025 at 04:41:38 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/vimentin-monoclonal-mouse-antibody-vim-3736/>