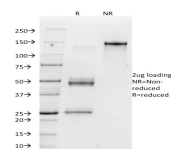


Vitronectin Receptor / CD51 / CD61 Monoclonal Mouse Antibody (23C6)



Product Description

ITAGV encodes integrin alpha chain V. Integrins are heterodimeric integral membrane proteins composed of an alpha chain and a beta chain. The I-domain containing integrin alpha V undergoes post-translational cleavage to yield disulfide-linked heavy and light chains, that combine with multiple integrin beta chains to form different integrins. Among the known associating beta chains (beta chains 1, 3, 5, 6, and 8; 'ITGB1', 'ITGB3', 'ITGB5', 'ITGB6', and 'ITGB8'), each can interact with extracellular matrix ligands; the alpha V beta 3 integrin, perhaps the most studied of these, is referred to as the Vitronectin Receptor (VNR). In addition to adhesion, many integrins are known to facilitate signal transduction. Alternative splicing results in multiple transcript variants. Reacts with Integrin Alpha V Beta 3. 23C6 may be useful for bone resorption modulation, osteoclast identification, receptor purification, malignant melanoma identification and treatment.

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 1471, Anti-CD51|CD61|Vitronectin Receptor (23C6)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#1471
Antibody reactivity (target)	CD51, CD61, Vitronectin Receptor
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	23C6
Isotype	IgG1, kappa
Molecular weight of antigen	150-160 kDa under non-reducing conditions. Under reducing conditions, it is cleaved into 2 bands of about 125-130 kDa +20-25 kDa.
Synonyms	CD51; DKFZp686A08142; integrin alpha V beta 3; Integrin alpha-5 Integrin alpha-V light chain; integrin alphaVbeta3; integrin; alpha V (vitronectin receptor; alpha polypeptide; antigen CD51); ITGAV; Msk8; VNRA; VTNR
Human gene symbol	ITGAV
Entrez gene ID	3685
SwissProt	P06756
Unigene	436873
Immunogen	Osteoclasts from osteoclastomas
Antibody target cellular localization	Plasma membrane
Species reactivity	Chicken, Human
Expected antibody applications	IHC (frozen) (published for clone), Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
Antibody application notes	ELISA: For coating, order antibody without BSA; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Squamous cells, A375M melanoma cells. A549 cells.
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
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.
Tumor expression	Melanoma

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.

References

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

1. J Cell Biol (1989) 109: 1817-1826. (IP, IF)
2. Exp Cell Res (1991) 195:368-375. (IP, functional studies, IF)
3. J Clin Invest (1992) 90: 1513-1522. (functional studies).
4. Am J Pathol (1997) 150(4): 1383-1393. (IHC, frozen)
5. Br J Cancer (1998) 77(4): 530-536. (Flow)

This datasheet was generated on June 17, 2026 at 10:09:31 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/vitronectin-receptor-cd51-cd61-monooclonal-mouse-antibody-23c6/>

