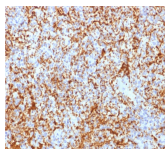


CD61 / Integrin Beta-3 / Platelet-Glycoprotein-IIIa Recombinant Monoclonal Rabbit Antibody (ITGB3/2166R)



Product Description

Reacts with human integrin beta3 (GPIIIa, vitronectin receptor beta chain). It associates with the IIb-chain (CD41) to form the GpIIb/GpIIIa complex (CD41/CD61). The CD41/CD61 complex appears early in megakaryocyte maturation. The activated CD41/CD61 complex is a receptor for von Willebrand factor, soluble fibrinogen, fibronectin, vitronectin and thrombospondin. It plays a central role in platelet activation and aggregation. The CD51/CD61 is implicated in tumor metastasis and adenoviral infection. The antibody detects platelets in smears of blood and bone marrow, as well as megakaryocytes in frozen sections and cell smears. The antibody is useful for classification of megakaryoblastic leukemia.

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 2166, Anti-CD61|Integrin 3|Platelet-Glycoprotein-IIIa (ITGB3/2166R)

Product attributes

Antibody number	#2166
Antibody reactivity (target)	CD61, Integrin 3, Platelet-Glycoprotein-IIIa
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	ITGB3/2166R
Isotype	IgG
Molecular weight	105 kDa & 90 kDa
Synonyms	BDPLT2; GP3A; GPIIIa; GT; HPA 1; HPA 4; Integrin beta-3; ITGB3; NAIT; Platelet fibrinogen receptor beta subunit; Platelet glycoprotein IIIa; PTP
Human gene symbol	ITGB3
Entrez gene ID	3690
SwissProt	P05106
Unigene	218040
Immunogen	Recombinant human ITGB3 protein fragment (aa 385-490) (exact sequence is proprietary)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	U937, KG1a, HEL cells. Spleen
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.
Cell/tissue expression	Platelets, Megakaryocytes
Antibody research areas	Cell adhesion
Tumor expression	Leukemia/lymphoma

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 December 11, 2025 at 02:39:28 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd61-integrin-beta-3-platelet-glycoprotein-iii-a-recombinant-monoclonal-rabbit-antibody-itgb3-2166r/>