

CD61 / Integrin Beta-3 / Platelet-Glycoprotein-Illa Monoclonal Mouse Antibody (25E11)

Product Description

Reacts with human integrin beta3 (GPIIIa, vitronectin receptor beta chain). The protein detectable is a complex of CD41 and CD61. The apparent molecular weight of the GPIIIa by SDS-PAGE is 105 kDa reduced and 90 kDa unreduced. Ligands are fibronectin, fibrinogen, von Willebrand factor, vitronectin and thrombospondin. Residues 237-248 of GPIIIa or CD61 are critical in adhesive protein binding.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL. Catalog number key for antibody number 1048, Anti-CD61 (25E11)

Product attributes

Product attributes				
Antibody number	#1048			
Antibody reactivity (target)	CD61			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	25E11			
Isotype	IgM, kappa			
Molecular weight	105 kDa & 90 kDa			
Synonyms	BDPLT2; GP3A; GPIlla; GT; HPA 1; HPA 4; Integrin beta-3; ITGB3; NAIT; Platelet fibrinogen receptor beta subunit; Platelet glycoprotein Illa; PTP			
Human gene symbol	ITGB3			
Entrez gene ID	3690			
SwissProt	P05106			
Unigene	218040			
Immunogen	Normal blood mononuclear cells activated by mixed lymphocyte reaction			
Antibody target cellular localization	Plasma membrane			
	Human			
Species reactivity	Human			
Species reactivity Antibody application notes	Human For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry			
, ,	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/ml. for most applications, or 1 ug/million			
Antibody application notes	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or			
Antibody application notes Positive control	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils.			
Antibody application notes Positive control Shipping condition	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils. Room temperature Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35			
Antibody application notes Positive control Shipping condition Storage Conditions	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils. Room temperature Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils. Room temperature Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils. Room temperature Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% aside, Purified: BSA-free: 1 mg/mL in			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils. Room temperature Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 1 mg/mL in PBS without azide			

Call us: 800-304-5357 Email: techsupport@biotium.com

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

McMichael AJ et al. (eds) Leukocyte Typing III, Oxford University Press, Oxford, 1987. | Schlossman S. et al. (eds) Leukocyte Typing V, Oxford University Press, Oxford, 1995

This datasheet was generated on October 30, 2025 at 04:07:35 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd61-25e11/