



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

CD61 / Integrin Beta-3 / Platelet-Glycoprotein-IIIa 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.

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

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; 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	Normal blood mononuclear cells activated by mixed lymphocyte reaction
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
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
Positive control	U937, KG1a, HEL cells, and human platelets in lymph nodes or tonsils.
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/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
Cell/tissue expression	Platelets, Megakaryocytes
Antibody research areas	Cell adhesion

This datasheet was generated on July 20, 2024 at 08:26:07 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-cd61-25e11/>