CD57 / B3GAT1 Monoclonal Mouse Antibody (NK-1 + HNK-1)



Product Description

Anti-CD57 marks a subset of lymphocytes known as natural killer (NK) cells. Follicular center cell lymphomas often contain many NK cells within the neoplastic follicles. Anti-CD57 also stains neuroendocrine cells and their derived tumors, including carcinoid tumor and medulloblastoma. Anti-CD57 can also be useful in separating type B3 thymoma from thymic carcinoma when combined with a panel that includes antibodies against GLUT1, CD5, and CEA.

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 0695, Anti-B3GAT1|CD57 (NK-1 HNK-1)

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

Product attributes

Antibody number	#0695		
Antibody reactivity (target)	B3GAT1, CD57		
Antibody type	Primary		
Host species	Mouse		
Clonality	use noclonal -1 + HNK-1 M 10 kDa Slucuronyltransferase 1; B3GAT1; lactosylgalactosylxylosylprotein beta-Glucuronosyltransferase 1; GLCATP; GlcUAT-P; lcuronosyltransferase P; UDP GlcUA Glycoprotein beta 1, 3 lcuronyltransferase GAT1 087 P2W7 1050 man peripheral blood mononuclear cells (HNK-1 & NK-1) lisma membrane C (FFPE) (verified) man pher concentration may be required for direct detection using mary antibody conjugates than for indirect detection with, bondary antibody, Immunofluorescence: 0.5-1 ug/ml., Does		
Clone	NK-1 + HNK-1		
Isotype	IgM		
Molecular weight	~110 kDa		
Synonyms	3-Glucuronyltransferase 1; B3GAT1; Galactosylgalactosylxylosylprotein 3-beta-Glucuronosyltransferase 1; GLCATP; GlcUAT-P; Glucuronosyltransferase P; UDP GlcUA Glycoprotein beta 1, 3 Glucuronyltransferase		
Human gene symbol	B3GAT1		
Entrez gene ID	27087		
SwissProt	Q9P2W7		
Unigene	381050		
Immunogen	Human peripheral blood mononuclear cells (HNK-1 & NK-1)		
Antibody target cellular localization	Plasma membrane		
Verified antibody applications	IHC (FFPE) (verified)		
Species reactivity	Human		
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mt_, Does not react with rat; others not known, Immunohistology formalin-fixed 2-4 ug/mt_, Staining of formalin-fixed dissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user		
Positive control	Lymph node or Tonsil		
Shipping condition	Room temperature		
Storage Conditions	Store at 2 to 8 $^{\circ}\text{C}$, Note: store BSA-free antibodies at -10 to -35 $^{\circ}\text{C}$		
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		
Antibody research areas	Cancer, Immunology		
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 October 20, 2025 at 05:10:24 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd57-b3gat1-nk-1-hnk-1/