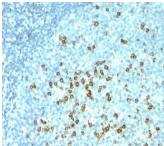


CD57 / B3GAT1 Monoclonal Mouse Antibody (NK-1 + HNK-1)



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.

Product attributes	
Antibody number	#0695
Antibody reactivity (target)	B3GAT1, CD57
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
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/mL, Does not react with rat; others not known, Immunohistology formalin-fixed 2-4 ug/mL, Staining of formalin-fixed tissues 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 °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
Antibody research areas	Cancer, Immunology