

BD Biosciences Mouse CD Marker Chart



bdbiosciences.com/go/mousedmarkers

CD	Alternative Name	Ligands & Associated Molecules	T Cell	B Cell	Dendritic Cell	Macrophage	Granulocyte	Platelet	Endothelial Cell	Epithelial Cell
CD16	CD1, CD2, Ly-38	Lipid, Glycolipid Ag	+	+	+	+	+	+	+	+
CD2	UFA-2, Ly-37, Ly-37	CD48, CD58, CD59, CD15	+	+	+	+	+	+	+	+
CD3	CD3 γ , CD3 δ , TCR γ	TCR complex	+	+	+	+	+	+	+	+
CD3a	CD3 α , CD3 δ , TCR α	TCR complex	+	+	+	+	+	+	+	+
CD3b	CD3 β , CD3 δ , TCR β	TCR complex	+	+	+	+	+	+	+	+
CD4	CD4, Ly-4	MHC class II, HV gp120, IL-16	+	+	+	+	+	+	+	+
CD5	Ly-1, Ly-11, Ly-12, Ly-A	CD72	+	+	+	+	+	+	+	+
CD5.1	Ly-1.1	CD72	+	+	+	+	+	+	+	+
CD6	T12	CD166 (ALCAM), BA11	+	+	+	+	+	+	+	+
CD7	gp40		+	+	+	+	+	+	+	+
CD8a	Ly-2, Ly-2.2, Ly-8, Ly-35	MHC class I	+	+	+	+	+	+	+	+
CD8b	Ly-3, Ly-3.2, Ly-4C, CD8b1	MHC class I	+	+	+	+	+	+	+	+
CD8b.2	Ly-3.2, Ly-8.2	MHC class I	+	+	+	+	+	+	+	+
CD9	Tapanzi		+	+	+	+	+	+	+	+
CD10	CALLA, MME, NRP	Proteoglycans	+	+	+	+	+	+	+	+
CD10a	Ly-15, Ly-21, Integrin α L	CD54, CD102, CD50	+	+	+	+	+	+	+	+
CD10b	Integrin α L, Ly-40, CD3, CD1A, MHC I	CD54, ICAM-1, Fibronectin	+	+	+	+	+	+	+	+
CD11c	ITGA6 (Integrin α X), CD4 (antigen receptor α), ICAM-1 receptor, Ly-6G, Ly-6S, Ly-6F, CD11c	ICAM-1, Fibronectin, ICAM-1	+	+	+	+	+	+	+	+
CD117	Aminopeptidase N, gp150	L-Asparaginase	+	+	+	+	+	+	+	+
CD14	M2, I-PS Receptor	LPS/PS complex, TLR2, TLR4	+	+	+	+	+	+	+	+
CD15	SSEA-1, F4/20, Lewis x	CD62	+	+	+	+	+	+	+	+
CD16	Fc γ RIII, Fc γ RIIIa, Ly-17, FCGR3, FCGR3	IGFC	+	+	+	+	+	+	+	+
CD18	Integrin β 2	CD11a, b, c	+	+	+	+	+	+	+	+
CD19	B4	CD19, CD81	+	+	+	+	+	+	+	+
CD20	Ly-44, B1		+	+	+	+	+	+	+	+
CD21	CD21/22	C8E, EBV, CD23, CD19, CD81	+	+	+	+	+	+	+	+
CD22	Ly-8.2, Sigslec-2	N-Glycolyl neuraminic acid	+	+	+	+	+	+	+	+
CD23	Fc γ RII, Ly-42	IgE, CD21, CD11a, CD11c	+	+	+	+	+	+	+	+
CD24	Heat Stable Antigen, Ly-52, Hectadlin	CD69 (P-selectin)	+	+	+	+	+	+	+	+
CD25	Ly-43, IL-2 Receptor α chain, p25	IL-2 Receptor α	+	+	+	+	+	+	+	+
CD26	Dipeptidyl peptidase, DPP IV, TMSM	Peptidases	+	+	+	+	+	+	+	+
CD27	T14, v152, v167, T155		+	+	+	+	+	+	+	+
CD28	T1504 antigen or T1544	CD80, B7-1, CD86, B7-2	+	+	+	+	+	+	+	+
CD29	Integrin β 1, VLA5, gp130	VCAM-1, MAdCAM-1, ECM	+	+	+	+	+	+	+	+
CD30	Ki-1	CD153	+	+	+	+	+	+	+	+
CD31	PECAM-1, gPIIb, mAbCAM, platelet endothelial cell adhesion molecule, PECAM-1	CD31, Vitronectin receptor	+	+	+	+	+	+	+	+
CD32	Fc γ RII, Ly-17, Ly-122, Fc γ receptor 2, Low affinity immunoglobulin γ 2c receptor	IgG	+	+	+	+	+	+	+	+
CD33	gp87, M2E1C, Sialic acid binding by like lectin II, myeloid cell surface antigen CD33	Sialylated glycoproteins, Sugar chains containing sialic acid, sialic acid 2-6-linked sialic acid	+	+	+	+	+	+	+	+
CD34	Mucosialin	CD46 (L-selectin)	+	+	+	+	+	+	+	+
CD35	CR1, CR3 receptor	CR1, CR3, CR3a	+	+	+	+	+	+	+	+
CD36	Scavenger receptors (SR), SRB, SRB3	Oxidized LDL, Thrombospondin, Collagen	+	+	+	+	+	+	+	+
CD37	gp130-42, Leukocyte antigen CD37, Teneurin-2b, TSPAN5	CD35, CD81, CD82, MHC class II	+	+	+	+	+	+	+	+
CD38	ADP-ribosyl cyclase, T10, Cyclic ADP-ribose hydrolase 1	CD31, Hyaluronic acid, CD37-like complex, CD16, MHC class II	+	+	+	+	+	+	+	+
CD39	NTF50a, E. Faecalibacterium prausnitzii, diacylglycerol transferase 1 (DAGAT1), ATP1B2, hydroxymethylglutaryl-CoA lyase, HMG-CoA lyase	ATP, ADP	+	+	+	+	+	+	+	+
CD40	gp39 receptor, gp39, MGC0013, TNFRSF3, Tumor necrosis factor receptor superfamily member 5	ATP, ADP	+	+	+	+	+	+	+	+
CD41	GPIIb, Integrin α IIb, Platelet membrane glycoprotein IIb, ITGA2b, Integrin α IIb, human Platelet Antigen-2 (HPA-2)	Fibronectin, Fibrinogen, von Willebrand factor, Thrombospondin	+	+	+	+	+	+	+	+
CD42a	GPIIb, Platelet glycoprotein IIb	von Willebrand factor	+	+	+	+	+	+	+	+
CD42b	GPIIb, Platelet glycoprotein IIb	von Willebrand factor	+	+	+	+	+	+	+	+
CD42c	GPIIb, Platelet glycoprotein IIb	von Willebrand factor	+	+	+	+	+	+	+	+
CD43	Ly-48, Sialoprotein, Leukosialin, Galactosyltransferase, SPN	CD43a	+	+	+	+	+	+	+	+
CD44	Ly-24, E-Cadherin, H-CAM, Pgp-1, Phagocyte glycoprotein I, Extracellular matrix receptor, Ly-24, lymphocyte homing/adhesion receptor, Myxomatous receptor	Fibronectin, Collagen, Fibronectin, Laminin, Osteopontin	+	+	+	+	+	+	+	+
CD45	Leukocyte Common Antigen (LCA)	CD135, CD11c, CD22, CD4, CD45AP, p56lck, p59fyn, Src kinases	+	+	+	+	+	+	+	+
CD45.1	Ly-5.1, Ly-5, P-ITRC		+	+	+	+	+	+	+	+
CD45.2	Ly-5.2, Ly-5b, P-ITRC		+	+	+	+	+	+	+	+
CD45R	B220, Ly-5.1, Ly-4, T200, Protein tyrosine phosphatase receptor type C (PTP-RC)		+	+	+	+	+	+	+	+
CD45RB			+	+	+	+	+	+	+	+
CD45RC	Protein tyrosine phosphatase, receptor type C (PTP-RC)		+	+	+	+	+	+	+	+
CD45RO	Membrane Cell Surface Protein (MCP), T-lymphocyte leukocyte common antigen, TR42.10	CR3, CR4, Molecules virus	+	+	+	+	+	+	+	+
CD46	Integrin-associated protein (IAP), OAK3, Neuropilin, MEB, gp42	SERP (CD172), CD81, Thrombospondin	+	+	+	+	+	+	+	+
CD48	Ban-1, HLA-8, BCM-1, DX-45, MEM-102	CD2, ICAM-1, CD229, CD244	+	+	+	+	+	+	+	+
CD49a	VLA-1a, Integrin α 1	Collagen, Laminin	+	+	+	+	+	+	+	+
CD49b	VLA-2a, Integrin α 2, gp7a	Collagen, Laminin, MMP-1	+	+	+	+	+	+	+	+
CD49c	VLA-3a, Integrin α 3, GARP3, Galactosyltransferase B3, M6E18, Very Common Antigen-2 (VCA-2)	Fibronectin, laminin, collagen	+	+	+	+	+	+	+	+
CD49d	VLA-4a, Integrin α 4	CD106 (VCAM1), MAdCAM, Fibronectin, Plectin	+	+	+	+	+	+	+	+
CD49e	VLA-5a, Integrin α 5, Fibronectin receptor	Fibronectin, Ixvactin, Fibrinogen	+	+	+	+	+	+	+	+
CD49f	VLA-6a, Integrin α 6, gp7	Laminin, Ixvactin	+	+	+	+	+	+	+	+
CD51	Integrin α 5, VLA-6a, Vitronectin α 5	Vitronectin, Fibronectin, Fibrinogen, thrombospondin, von Willebrand factor, CD31	+	+	+	+	+	+	+	+
CD52	CAMPATH-1, HEA, Epitaxial secondary protein B2, CLS1, MB7, B7	VCAM-1, Integrins	+	+	+	+	+	+	+	+
CD53	MOXA4, TSPAN25, TSPAN25	VCAM-1, Integrins	+	+	+	+	+	+	+	+
CD54	ICAM-1, MAdCAM-2	ICAM-1, MAdCAM-1, CD31, CD11a/CD11b, Rhesus, CD227	+	+	+	+	+	+	+	+
CD55	Decoy and signaling factor (DAG), complement-glycosylphosphatidylinositol, Complement blood group, Sd-50, DAF, GPI-DAF	CR1, CR3, CD91, CD225, CD315, CD316	+	+	+	+	+	+	+	+
CD56	Ly-19, N19, Neural Cell Adhesion Molecule (NCAM)	NCAM-1, Heparin sulfate	+	+	+	+	+	+	+	+
CD57	T154g, N19, sialoside, MACS, MRL-1, P18	CR3, CR4, ICAM-1	+	+	+	+	+	+	+	+
CD61	GPIIb, Integrin β 3	Fibrinogen, Fibrin, TGF β , TSP1 and CR1	+	+	+	+	+	+	+	+
CD62E	E-selectin, ELAM-1, LECAM-2	Sialyl Lewis x, CLA, CD62	+	+	+	+	+	+	+	+
CD62L	L-selectin, LECAM-1, Lfne, Ly-22, Ly-22, Ly-22, Ly-22, Ly-22, Ly-22, Ly-22	CD64, MHC class II	+	+	+	+	+	+	+	+
CD62P	P-selectin, GMP-140, PADGEM	CD62L, CD24	+	+	+	+	+	+	+	+
CD63	LAMP, M6A1, gp55, M6A, LAMP-3, M6A91, OMAH8, TSPAN10, granule/lysosomal membrane 1 antigen	VCAM-1, VLA-6, CD81, CD8, PR4 kinase, CD117, CD82	+	+	+	+	+	+	+	+
CD64	Fc γ RI, Fc γ receptor 1, High affinity immunoglobulin γ 1c receptor 1 (Fc γ RIa)	IgG	+	+	+	+	+	+	+	+
CD66a	B22, CD66a		+	+	+	+	+	+	+	+
CD66b	CD66b, CD66b		+	+	+	+	+	+	+	+
CD68	Macrosialin (lysosomal glycoprotein 110)	LDL	+	+	+	+	+	+	+	+
CD69	Very Early Adhesion Antigen		+	+	+	+	+	+	+	+
CD70	CD27 Ligand	CD27	+	+	+	+	+	+	+	+
CD71	Transferrin Receptor	Transferrin	+	+	+	+	+	+	+	+
CD72	gp-2, Ly-119	CD5, CD100	+	+	+	+	+	+	+	+
CD73	HEV, Ecto-5'-nucleotidase	MMP	+	+	+	+	+	+	+	+
CD74	NK associated invariant chain (i)	MHC class II	+	+	+	+	+	+	+	+
CD77	PK blood group antigen (B), CTR, Gb3, α -1,4-Galactosyltransferase, Gb3 synthase, Agp40		+	+	+	+	+	+	+	+
CD79a	Ig α , μ 1, Ly-54	Ig α , CD5, CD19, CD279	+	+	+	+	+	+	+	+
CD79b	Ig β , B29	Ig β , CD5, CD19, CD279	+	+	+	+	+	+	+	+
CD80	B7-1, B7-1, Ly-33	CD28, CD152	+	+	+	+	+	+	+	+
CD81	TAPA-1	CD95, CD137, CD225, CD315, CD316	+	+	+	+	+	+	+	+
CD82	CD3 Ag, KH1	MHC molecules, CD4, CD8, CD20, CD37, CD81, Integrins	+	+	+	+	+	+	+	+
CD83	H815		+	+	+	+	+	+	+	+
CD84	G86	IL8	+	+	+	+	+	+	+	+
CD86	B7-2, B7.2, Ly-38	CD28, CD152	+	+	+	+	+	+	+	+
CD87	vLA receptor	vLA, Vitronectin	+	+	+	+	+	+	+	+
CD88	C5a ligand, C5aR	C5a	+	+	+	+	+	+	+	+
CD90	Thy-1, α , T25		+	+	+	+	+	+	+	+
CD90.1	Thy-1.2, α MR		+	+	+	+	+	+	+	+
CD90.2	Thy-1.2, α CH		+	+	+	+	+	+	+	+
CD93	LRF, A2M, AD185f2, α 2-Macroglobulin receptor	LDL, LRPAP1, α 2M	+	+	+	+	+	+	+	+
CD94	4A4.1, C1qR	CD11	+	+	+	+	+	+	+	+
CD94	KP43, hml1	IgD, IgM	+	+	+	+	+	+	+	+
CD95	Fas, APO-1	CD178 (Fas ligand)	+	+	+	+	+	+	+	+
CD96	Tacile	CD135, Nectin-1	+	+	+	+	+	+	+	+
CD97	TM13b1, TMS5	CD55	+	+	+	+	+	+	+	+
CD98	4F2, Ly-18, R388	CD28, CD147, Topoisomerase, Actin	+	+	+	+	+	+	+	+

CD	Alternative Name	Ligands & Associated Molecules	T Cell	B Cell	Dendritic Cell	Macrophage	Granulocyte	Platelet	Endothelial Cell	Epithelial Cell
CD99	Paired immunoglobin-like type 2 receptor ligand	IL18R	+	+	+	+	+	+	+	+
CD100	Semaphorin H, col-4A, Semaphorin 4D, Semu 4D	CD72, Plexin-B1, CD45	+	+	+	+	+	+	+	+
CD101	IG2F, V7		+	+	+	+	+	+	+	+
CD102	ICAM-2, 40	ICAM-1, Mac-1	+	+	+	+	+	+	+	+
CD103	Integrin α L	E-Cadherin	+	+	+	+	+	+	+	+
CD104	Integrin β 4	Laminin, Plectin	+	+	+	+	+	+	+	+
CD105	Endoglin	TGF β	+	+	+	+	+	+	+	+
CD106	VCAM-1	VLA-4	+	+	+	+	+	+	+	+
CD107a	LAMP-1	Collagen, Laminin, Fibronectin	+	+	+	+	+	+	+	+
CD107b	LAMP-2, LAMP-2B, LAMP-8		+	+	+	+	+	+	+	+
CD108	Sema7a	Tyrosine kinases	+	+	+	+	+	+	+	+
CD110	Thrombospondin, Receptor c1orf	JAK2, Thrombospondin	+	+	+	+	+	+	+	+
CD111	PR1, Nectin-1, CD111, CD111, ED4, HgR, HVC, MGC142031, MGC142037, OC7, PR8, PR9, PR11, Sx-12	α -Herpesvirus, Nectin-3, Adelin	+	+	+	+	+	+	+	+
CD112	PR2, Nectin-2, PR2-2, HV8B, MPH, Psc, CD112, A232526, A987993, Receptor 2, PR2	CD226, PR8, α Adelin	+	+	+	+	+	+	+	+
CD113	PVR1, Nectin-3	Nectin-1, Nectin-2, PVR	+	+	+	+	+	+	+	+
CD114	G-CSF Receptor, G-CSFR, G-CSFR, Granulocyte colony-stimulating factor receptor, G-CSFR	G-CSF, JAK2, JAK2	+	+	+	+	+	+	+	+
CD115	M-CSF Receptor, CSF-1R, c-fms, Fms-2	M-CSF	+	+	+	+	+	+	+	+
CD116	GM-CSF Receptor α chain	GM-CSF	+	+	+	+	+	+	+	+
CD117	c-kit, Steel factor receptor, Dominant white spotting	Ligand Receptor, Stem cell factor, or mast-cell growth factor	+	+	+	+	+	+	+	+
CD118	IFN- γ Receptor, Type 1 IFN- γ , IFN- γ Receptor	IFN- γ , IFN- β	+	+	+	+	+	+	+	+
CD119	IFN- γ Receptor α chain	IFN- γ	+	+	+	+	+	+	+	+
CD120a	TNFR1, TNF-R55	TNF (Lymphostatin A (TNF- α))	+	+	+	+	+	+	+	+
CD120b	TNFR2, TNF-R75	TNF (Lymphostatin A (TNF- α))	+	+	+	+	+	+	+	+
CD121a	IL-1 Receptor Type 1	IL-1 α , IL-1 β	+	+	+	+	+	+	+	+
CD121b	IL-1 Receptor Type 2	IL-1 α , IL-1 β	+	+	+	+	+	+	+	+
CD122	IL-2 and IL-15 Receptor β chain	IL-2, IL-15	+	+	+	+	+	+	+	+
CD123	IL-3 Receptor α chain	IL-3	+	+</						