

BD Biosciences Human CD Marker Chart

Reviewed by HLDA¹
bdbiosciences.com/cdmarkers



23-12999-01

CD	Alternative Name	Ligands & Associated Molecules	T Cell	B Cell	Dendritic Cell	MK Cell	Stem Cell/Progenitor	Macrophage/Monocyte	Platelet	Granulocyte	Erythrocyte	Endothelial Cell	Epithelial Cell
CD1	CD1		-	-	-	-	-	-	-	-	-	-	-
CD2	CD2		+	+	+	+	+	+	+	+	+	+	+
CD3	CD3		+	+	+	+	+	+	+	+	+	+	+
CD4	CD4		+	+	+	+	+	+	+	+	+	+	+
CD5	CD5		+	+	+	+	+	+	+	+	+	+	+
CD6	CD6		+	+	+	+	+	+	+	+	+	+	+
CD7	CD7		+	+	+	+	+	+	+	+	+	+	+
CD8	CD8		+	+	+	+	+	+	+	+	+	+	+
CD9	CD9		+	+	+	+	+	+	+	+	+	+	+
CD10	CD10		+	+	+	+	+	+	+	+	+	+	+
CD11	CD11		+	+	+	+	+	+	+	+	+	+	+
CD12	CD12		+	+	+	+	+	+	+	+	+	+	+
CD13	CD13		+	+	+	+	+	+	+	+	+	+	+
CD14	CD14		+	+	+	+	+	+	+	+	+	+	+
CD15	CD15		+	+	+	+	+	+	+	+	+	+	+
CD16	CD16		+	+	+	+	+	+	+	+	+	+	+
CD17	CD17		+	+	+	+	+	+	+	+	+	+	+
CD18	CD18		+	+	+	+	+	+	+	+	+	+	+
CD19	CD19		+	+	+	+	+	+	+	+	+	+	+
CD20	CD20		+	+	+	+	+	+	+	+	+	+	+
CD21	CD21		+	+	+	+	+	+	+	+	+	+	+
CD22	CD22		+	+	+	+	+	+	+	+	+	+	+
CD23	CD23		+	+	+	+	+	+	+	+	+	+	+
CD24	CD24		+	+	+	+	+	+	+	+	+	+	+
CD25	CD25		+	+	+	+	+	+	+	+	+	+	+
CD26	CD26		+	+	+	+	+	+	+	+	+	+	+
CD27	CD27		+	+	+	+	+	+	+	+	+	+	+
CD28	CD28		+	+	+	+	+	+	+	+	+	+	+
CD29	CD29		+	+	+	+	+	+	+	+	+	+	+
CD30	CD30		+	+	+	+	+	+	+	+	+	+	+
CD31	CD31		+	+	+	+	+	+	+	+	+	+	+
CD32	CD32		+	+	+	+	+	+	+	+	+	+	+
CD33	CD33		+	+	+	+	+	+	+	+	+	+	+
CD34	CD34		+	+	+	+	+	+	+	+	+	+	+
CD35	CD35		+	+	+	+	+	+	+	+	+	+	+
CD36	CD36		+	+	+	+	+	+	+	+	+	+	+
CD37	CD37		+	+	+	+	+	+	+	+	+	+	+
CD38	CD38		+	+	+	+	+	+	+	+	+	+	+
CD39	CD39		+	+	+	+	+	+	+	+	+	+	+
CD40	CD40		+	+	+	+	+	+	+	+	+	+	+
CD41	CD41		+	+	+	+	+	+	+	+	+	+	+
CD42	CD42		+	+	+	+	+	+	+	+	+	+	+
CD43	CD43		+	+	+	+	+	+	+	+	+	+	+
CD44	CD44		+	+	+	+	+	+	+	+	+	+	+
CD45	CD45		+	+	+	+	+	+	+	+	+	+	+
CD46	CD46		+	+	+	+	+	+	+	+	+	+	+
CD47	CD47		+	+	+	+	+	+	+	+	+	+	+
CD48	CD48		+	+	+	+	+	+	+	+	+	+	+
CD49	CD49		+	+	+	+	+	+	+	+	+	+	+
CD50	CD50		+	+	+	+	+	+	+	+	+	+	+
CD51	CD51		+	+	+	+	+	+	+	+	+	+	+
CD52	CD52		+	+	+	+	+	+	+	+	+	+	+
CD53	CD53		+	+	+	+	+	+	+	+	+	+	+
CD54	CD54		+	+	+	+	+	+	+	+	+	+	+
CD55	CD55		+	+	+	+	+	+	+	+	+	+	+
CD56	CD56		+	+	+	+	+	+	+	+	+	+	+
CD57	CD57		+	+	+	+	+	+	+	+	+	+	+
CD58	CD58		+	+	+	+	+	+	+	+	+	+	+
CD59	CD59		+	+	+	+	+	+	+	+	+	+	+
CD60	CD60		+	+	+	+	+	+	+	+	+	+	+
CD61	CD61		+	+	+	+	+	+	+	+	+	+	+
CD62	CD62		+	+	+	+	+	+	+	+	+	+	+
CD63	CD63		+	+	+	+	+	+	+	+	+	+	+
CD64	CD64		+	+	+	+	+	+	+	+	+	+	+
CD65	CD65		+	+	+	+	+	+	+	+	+	+	+
CD66	CD66		+	+	+	+	+	+	+	+	+	+	+
CD67	CD67		+	+	+	+	+	+	+	+	+	+	+
CD68	CD68		+	+	+	+	+	+	+	+	+	+	+
CD69	CD69		+	+	+	+	+	+	+	+	+	+	+
CD70	CD70		+	+	+	+	+	+	+	+	+	+	+
CD71	CD71		+	+	+	+	+	+	+	+	+	+	+
CD72	CD72		+	+	+	+	+	+	+	+	+	+	+
CD73	CD73		+	+	+	+	+	+	+	+	+	+	+
CD74	CD74		+	+	+	+	+	+	+	+	+	+	+
CD75	CD75		+	+	+	+	+	+	+	+	+	+	+
CD76	CD76		+	+	+	+	+	+	+	+	+	+	+
CD77	CD77		+	+	+	+	+	+	+	+	+	+	+
CD78	CD78		+	+	+	+	+	+	+	+	+	+	+
CD79	CD79		+	+	+	+	+	+	+	+	+	+	+
CD80	CD80		+	+	+	+	+	+	+	+	+	+	+
CD81	CD81		+	+	+	+	+	+	+	+	+	+	+
CD82	CD82		+	+	+	+	+	+	+	+	+	+	+
CD83	CD83		+	+	+	+	+	+	+	+	+	+	+
CD84	CD84		+	+	+	+	+	+	+	+	+	+	+
CD85	CD85		+	+	+	+	+	+	+	+	+	+	+
CD86	CD86		+	+	+	+	+	+	+	+	+	+	+
CD87	CD87		+	+	+	+	+	+	+	+	+	+	+
CD88	CD88		+	+	+	+	+	+	+	+	+	+	+
CD89	CD89		+	+	+	+	+	+	+	+	+	+	+
CD90	CD90		+	+	+	+	+	+	+	+	+	+	+
CD91	CD91		+	+	+	+	+	+	+	+	+	+	+
CD92	CD92		+	+	+	+	+	+	+	+	+	+	+
CD93	CD93		+	+	+	+	+	+	+	+	+	+	+
CD94	CD94		+	+	+	+	+	+	+	+	+	+	+
CD95	CD95		+	+	+	+	+	+	+	+	+	+	+
CD96	CD96		+	+	+	+	+	+	+	+	+	+	+
CD97	CD97		+	+	+	+	+	+	+	+	+	+	+
CD98	CD98		+	+	+	+	+	+	+	+	+	+	+
CD99	CD99		+	+	+	+	+	+	+	+	+	+	+
CD100	CD100		+	+	+	+	+	+	+	+	+	+	+
CD101	CD101		+	+	+	+	+	+	+	+	+	+	+
CD102	CD102		+	+	+	+	+	+	+	+	+	+	+
CD103	CD103		+	+	+	+	+	+	+	+	+	+	+
CD104	CD104		+	+	+	+	+	+	+	+	+	+	+
CD105	CD105		+	+	+	+	+	+	+	+	+	+	+
CD106	CD106		+	+	+	+	+	+	+	+	+	+	+
CD107	CD107		+	+	+	+	+	+	+	+	+	+	+
CD108	CD108		+	+	+	+	+	+	+	+	+	+	+
CD109	CD109		+	+	+	+	+	+	+	+	+	+	+
CD110	CD110		+	+	+	+	+	+	+	+	+	+	+
CD111	CD111		+	+	+	+	+	+	+	+	+	+	+
CD112	CD112		+	+	+	+	+	+	+	+	+	+	+
CD113	CD113		+	+	+	+	+	+	+	+	+	+	+
CD114	CD114		+	+	+	+	+	+	+	+	+	+	+
CD115	CD115		+	+	+	+	+	+	+	+	+	+	+
CD116	CD116		+	+	+	+	+	+	+	+	+	+	+
CD117	CD117		+	+	+	+	+	+	+	+	+	+	+
CD118	CD118		+	+	+	+	+	+	+	+	+	+	+
CD119	CD119		+	+	+	+	+	+	+	+	+	+	+
CD120	CD120		+	+	+	+	+	+	+	+	+	+	+
CD121	CD121		+	+	+	+	+	+	+	+	+	+	+
CD122	CD122		+	+	+	+	+	+	+	+	+	+	+
CD123	CD123		+	+	+	+	+	+	+	+	+	+	+
CD124	CD124		+	+	+	+	+	+	+	+	+	+	+
CD125	CD125		+	+	+	+	+	+	+	+	+	+	+
CD126	CD126		+	+	+	+	+	+	+	+	+	+	+
CD127	CD127		+	+	+	+	+	+	+	+	+	+	+
CD128	CD128		+	+	+	+	+	+	+	+	+	+	+
CD129	CD129		+	+	+	+	+	+	+	+	+	+	+
CD130	CD130		+	+	+	+	+	+	+	+	+	+	+
CD131	CD131		+	+	+	+	+	+	+	+	+	+	+
CD132	CD132		+	+	+	+	+	+	+	+	+	+	+
CD133	CD133		+	+	+	+	+	+	+	+	+	+	+
CD134	CD134		+	+	+	+	+	+	+	+	+	+	+
CD135	CD135		+	+	+	+	+	+	+	+	+	+	+
CD136	CD136		+	+	+	+	+	+	+	+	+	+	+
CD137	CD137		+	+	+	+	+	+	+	+	+	+	+
CD138	CD138		+	+	+	+	+	+	+	+	+	+	+
CD139	CD139		+	+	+	+	+	+	+	+	+	+	+
CD140	CD140		+	+	+	+	+	+	+	+	+	+	+
CD141	CD141		+	+	+	+	+	+	+	+	+	+	+
CD142	CD142		+	+	+	+	+	+	+	+	+	+	+
CD143	CD143		+	+	+	+	+	+	+	+	+	+	+
CD144	CD144		+	+	+	+	+	+	+	+	+	+	+
CD145	CD145		+	+	+	+	+	+	+	+	+	+	+
CD146	CD146		+	+	+	+	+	+	+	+	+	+	+
CD147	CD147		+	+	+	+	+	+	+	+	+	+	+
CD148	CD148		+	+	+	+	+	+	+	+	+	+	+
CD149	CD149		+	+	+	+	+	+	+	+	+	+	+
CD150	CD150		+	+	+	+	+	+	+	+	+	+	+
CD151	CD151		+	+	+	+	+	+	+	+	+	+	+
CD152	CD152		+	+	+</								