



CLINICAL SINGLE COLOR ANTIBODIES

US Version



ASR Single Color

Description	Clone	Laser Emission Maximum	Status	Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)					
				455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
				PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD1a	BL6	ASR	-	-	B25330	-	IM1942U	-	IM3610	-	-	IM3645	-	-	-	-	
CD1c	L161	ASR	-	-	-	-	-	-	-	B46036	-	-	-	-	-	-	
CD2	SFC13Pt2H9 (T11)	ASR	-	-	B36533	-	629776	-	-	-	-	-	-	-	-	-	
	39C1.5	ASR	B09685	-	IM0442U	-	IM0443U	B46023	IM2634U	-	A51074	A60794	-	-	B12111	B01681	
CD3	UCHT1	ASR	A93687	B00068	IM1281U	-	IM1282U	IM2705U	6607010 IM2635U	A66327	-	IM2467U	-	-	B10823	A66329	
CD4	SFC112T4D11 (T4)	ASR	-	-	-	-	-	6604727	-	-	-	-	-	-	-	-	
	13B8.2	ASR	A82789	A96417	IM0448U	-	IM0449U	-	IM2636U	B16491	-	IM2468U	-	-	B10824	A94685	
CD5	BL1a	ASR	A82790	-	IM0468U	-	IM0469U	A33096	IM2637U	A70203	A51075	A60790	-	-	A78835	A78836	
	CLB-T1/1	ASR	-	-	-	-	-	B55387	-	-	-	B55386	-	-	-	-	
CD6	2H46D3B	ASR	-	-	B16492	-	-	-	-	-	-	-	-	-	-	-	
CD7	3A1E-12H7	ASR	-	-	6603824 6603825	-	6603822 6603827	-	-	-	-	-	-	-	-	-	
	8H8.1	ASR	B06499	-	IM0585U	-	IM1429U	A70202	IM3613U	-	A46526	A97050	-	-	A70201	B16892	
CD8	SFC121Thy2D3 (T8)	ASR	-	-	-	-	-	6604728	6607011	A99019	-	A99023	-	-	-	-	
	B9.11	ASR	A82791	B00067	IM0451U	-	IM0452U	B08467	IM2638U	B21205	-	IM2469U	-	B76279	A66332	A94686	
CD8β	25T8.5H7	ASR	-	-	B42025	-	IM2217U	-	-	-	-	-	-	-	-	-	
CD9	ALB6	ASR	B09979	-	IM1755U	-	-	-	-	-	-	-	-	-	-	B13649	
CD10	J5	ASR	-	-	6604107 6602463	-	6604120 6603475	-	-	-	-	-	-	-	-	-	
	ALB1	ASR	-	-	IM2720U	-	IM1915U	IM3608U	IM2721U	B16490	A46527	IM3633	-	-	A86353	A89310	
	ALB2	ASR	-	-	IM0471U	-	-	-	-	-	-	-	-	-	-	-	
CD11a	25.3	ASR	-	-	IM0860U	-	IM1433U	-	-	-	-	-	-	-	-	-	
CD11b	94 (Mo1)	ASR	-	-	6602573	-	-	-	-	-	-	-	-	-	-	-	
	Bear1	ASR	B16891	-	IM0530U	-	IM2581U	-	IM3611	-	A54822	A87782	-	-	-	A97052	
CD11c	BU15	ASR	-	-	-	-	IM1760U	-	IM3707	B19719	A80249	B01680	-	-	B43304	-	
CD13	366 (MY7)	ASR	-	-	-	-	6602989	-	-	-	B19714	-	-	-	-	-	
	Immu103.44	ASR	-	-	-	-	-	A33097	IM2639U	A79389	A46528	A87783	-	-	-	-	
	SJ1D1	ASR	-	-	IM0778U	-	IM1427U	B24026	-	-	-	-	-	-	-	-	
CD14	322A-1	ASR	-	-	B36630	-	-	-	-	-	-	-	-	-	-	-	
	116	ASR	-	-	B36631	-	B36634	-	-	-	-	-	-	-	-	-	
	RM052	ASR	B00846	B01175	IM0645U	-	IM0650U	IM2707U	IM2640U	A70204	A22331	IM2580U	-	-	A99020	A86052	
CD15	80H5	ASR	A74775	B01176	IM1423U	-	IM1954U	-	IM2641U	-	-	-	-	-	-		
CD16	3G8	ASR	A82792	B00069	6604894 IM0814U	-	IM1238U	A33098	IM2642U	-	-	B00845	-	-	B20023	A66330	
CD16b	1D3	ASR	-	-	IM2353U	-	-	-	-	-	-	-	-	-	-	-	
CD18	7E4	ASR	-	-	IM1568U	-	IM1570U	-	-	-	-	-	-	-	-	-	
CD19	89B	ASR	-	-	B36534	-	B36536	-	-	-	-	-	-	-	-	-	
	HD237	ASR	-	-	-	-	-	6604551	-	-	-	-	-	-	-	-	
	J3-119	ASR	A86355	A96418	IM1284U	-	IM1285U	IM2708U	IM2643U	A66328	IM3628U	IM2470U	-	B76283	A78837	A78838	

 New

AF 647 : Alexa Fluor® 647

AF 700 : Alexa Fluor® 700

APC-A700 : APC-Alexa Fluor® 700

APC-A750 : APC-Alexa Fluor® 750

PB : Pacific Blue®

KrO : Krome Orange

ASR Single Color

Description	Clone	Laser Emission Maximum Status	Violet Laser (405 nm)		Blue Laser (488 nm)							Red Laser (640 nm)				
			455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
			PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD20	H299	ASR	-	-	B36626	-	B36625	-	-	-	-	-	-	-	-	
	FMC7	ASR	B08466	-	IM1364U	-	-	-	-	-	-	-	-	-	-	
	B9E9	ASR	A74777	-	IM1455U	-	IM1451U	IM3607U	IM2644U	B23134	IM3629U	A51076	-	B76287	B12112	A66331
CD21	BL13	ASR	B09982	-	IM0473U	-	A32536	-	-	-	-	-	-	-	-	
CD22	HD239	ASR	-	-	-	-	6604428	-	-	-	-	-	-	-	-	
	SJ10.1H11	ASR	B06498	-	IM0779U	-	IM1835U	B10245	IM3704	A80712	A74788	A60791	-	-	A89311	-
CD23	9P25	ASR	A93689	-	IM0529U	-	A33099	IM3609U	-	-	B12935	A69964	-	-	A70207	-
	HD50	ASR	-	-	-	-	6604426	-	-	-	-	-	-	-	-	
CD24	ALB9	ASR	-	-	-	-	IM1428U	B12699	IM2645	B23133	-	A87785	-	-	-	B10738
CD25	1HT44H3	ASR	-	-	-	-	6604422	-	-	-	-	-	-	-	-	-
	B1.49.9	ASR	A82944	-	IM0478U	-	IM0479U	-	IM2646U	A79386	A52882	B09684	-	-	A86356	B13979
CD26	4EL-1C7	ASR	-	-	-	-	IM2724	-	-	-	-	-	-	-	-	-
CD27	1A4CD27	ASR	B30635	-	-	-	IM2578	B26603	-	B21444	A54823	B09983	-	-	-	B12701
CD28	CD28.2	ASR	-	-	IM1236U	-	IM2071U	-	-	B24027	B23313	B10244	-	-	B12696	B08757
CD29	K20	ASR	-	-	IM0791U	-	-	-	-	-	-	-	-	-	-	-
CD30	HRS4	ASR	-	-	-	-	IM2033U	-	-	-	-	A87939	-	-	B42017	-
CD31	5.6E	ASR	B13035	-	IM1431U	-	-	-	-	-	-	-	-	-	-	-
CD32	2E1	ASR	-	B01177	-	-	IM1935	-	-	-	-	-	-	-	-	-
CD33	906	ASR	-	-	-	-	6604121 6603042	-	-	-	-	-	-	-	-	-
	D3HL60.251	ASR	-	-	IM1135U	-	IM1179U	-	IM2647U	A70198	A54824	IM2471U	-	-	-	A70200
CD34	581	ASR	-	-	IM1870U	-	IM1871U	IM2709U	IM2648U	-	A51077	IM2472U	-	-	A86354	A89309
	QBEnd10	ASR	-	-	-	-	IM1250U	-	IM2650U	-	-	-	-	-	-	-
	QBEnd10 + Immu133 + Immu409	ASR	-	-	-	-	IM1459U	-	-	-	-	-	-	-	-	-
CD36	FA6.152	ASR	B43302	-	IM0766U	-	-	-	-	-	-	A87786	-	-	B46022	-
CD37	BL14	ASR	-	-	IM0457	-	IM0458	-	-	-	-	-	-	-	-	-
CD38	T16	ASR	-	-	IM0775U	-	IM1832U	-	-	-	-	-	-	-	-	-
	LS198-4-3	ASR	B09683	-	-	-	IM2371U 6604928	A99022	IM2651U	A70205	A54189	A60792	-	-	B23489	A86049
CD39	BA54	ASR	-	-	-	-	B55384	-	-	B55385	-	-	-	-	-	-
CD40	MAB89	ASR	-	-	-	-	IM1936U	-	-	B20025	-	B30636	-	-	-	-
CD41	P2	ASR	-	-	IM0649U	-	IM1416U	-	-	-	-	B16894	-	-	-	-
	S222	ASR	-	-	IM1756U	-	-	-	-	-	-	-	-	-	-	-
CD42a	SZ1	ASR	-	-	IM1757U	-	-	-	-	-	-	-	-	-	-	-
CD42b	SZ2	ASR	-	-	IM0648U	-	IM1417U	-	-	-	-	B13980	-	-	-	-
CD43	DFT1	ASR	-	-	IM3264U	-	A32560	-	-	-	-	-	-	-	-	A89307
CD44	J.173	ASR	B37789	-	IM1219U	-	A32537	-	-	-	-	-	-	-	-	B30637

 New

AF 647 : Alexa Fluor* 647

AF 700 : Alexa Fluor* 700

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

ASR Single Color

Description	Clone	Laser Emission Maximum	Status	Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)					
				455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
				PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD45	ALB12	ASR	-	-	IM0647	-	-	-	-	-	-	-	-	-	-	-	
	Immu19.2	ASR	-	-	-	-	-	-	IM2652U	-	-	-	-	-	-	-	
	J33	ASR	A74765	A96416	IM0782U	-	IM2078U	IM2710U	IM2653U	A54139	IM3548U	IM2473U	A94246	-	A71117	A71119	
	KC56	ASR	-	-	6604104 6603838	-	6603839	-	-	-	-	-	-	-	-	-	
CD45RA	ALB11	ASR	-	-	IM0584U	-	IM1834U	-	-	-	-	-	-	-	-	-	
	2H4LDH11LDB9	ASR	A82946	-	-	-	-	IM2711U	-	-	B10821	B14807	-	B90396	-	A86050	
CD45R0	UCHL1	ASR	-	-	IM1247U	-	IM1307U	IM2712U	-	B30638	B13648	-	-	-	-	-	
CD48	J4.57	ASR	-	-	IM1837U	-	-	B06026	-	-	-	-	-	-	-	-	
CD49d	HP2/1	ASR	-	-	IM1404U	-	-	-	-	-	-	B01682	-	-	-	B16893	
CD49e	SAM1	ASR	-	-	IM1854U	-	-	-	-	-	-	-	-	-	-	-	
CD52	HI186	ASR	-	-	-	-	-	-	-	-	B68152	-	-	-	-	-	
CD54	84H10	ASR	-	-	IM0726U	-	IM1239U	-	-	-	-	-	-	-	-	-	
CD55	JS11KSC2.3	ASR	-	-	IM2725	-	IM2726	-	-	-	-	-	-	-	-	-	
CD56	N901	ASR	-	-	-	-	IM2073U 6603857 6603067	A82943	IM2654U	A79388	A51078	IM2474U	-	-	B10822	B46024	
CD57	NC1	ASR	A74779	-	IM0466U	-	-	-	-	-	-	-	-	-	-	-	
CD58	AICD58	ASR	-	-	IM1218U	-	IM1430	-	IM3702	-	-	IM3701	-	-	-	-	
CD59	P282E	ASR	-	-	IM3457U	-	-	-	-	-	-	-	-	-	-	-	
	MEM-43	ASR	-	-	-	-	B68140	-	-	-	-	-	-	-	-	-	
CD61	SZ21	ASR	-	-	IM1758U	-	IM3605	-	-	B21172	IM3716	-	-	-	-	-	
CD62L	DREG56	ASR	B30640	-	IM1231U	-	IM2214U	IM2713U	IM2655U	-	B30641	B30639	-	-	-	B26604	
CD62P	CLB-Thromb/6	ASR	-	-	-	-	IM1759U	-	-	-	-	-	-	-	-	-	
CD63	CLB-Gran/12	ASR	-	-	IM1165U	-	IM1914U	-	-	-	-	-	-	-	-	-	
CD64	22	ASR	B19718	-	IM1604U	-	IM3601U	A98434	IM3606U	-	B06025	-	-	-	-	A89308	
CD65	88H7	ASR	B08164	-	IM1654U	-	-	-	-	-	-	-	-	-	-	-	
CD66b	80H3	ASR	-	-	IM0531U	-	-	-	-	-	-	B15091	-	-	-	B08756	
CD66c	KOR-SA3544	ASR	-	-	IM2039U	-	IM2357U	-	-	-	-	-	-	-	-	-	
CD69	TP1.55.3	ASR	-	-	-	-	IM1943U	-	-	-	A80710	A80711	-	-	-	-	
CD71	YDJ1.2.2	ASR	-	-	IM0483U	-	IM2001U	-	-	-	-	-	-	-	A97051	A89313	
CD73	AD-2	ASR	-	-	-	-	B68176	-	-	-	-	-	-	-	-	-	
CD79a	HM47	ASR	-	-	-	-	IM2221U	-	IM3456U	B42018	-	A60793	-	-	-	-	
CD79b	CB3-1	ASR	-	-	-	-	IM1612U	-	-	B21445	-	A87784	-	-	-	-	
CD80	MAB104	ASR	-	-	IM1853U	-	IM1976U	-	-	-	B30644	B30642	-	-	-	B30643	
CD81	JS64	ASR	B19717	-	B25329	-	IM2579	-	-	-	-	A87789	-	-	-	-	
CD83	HB15a	ASR	-	-	IM2410U	-	IM2218U	-	IM3240U	-	-	-	-	-	-	-	
CD85j	HP-F1	ASR	-	-	-	-	A07408	-	-	-	-	B30645	-	-	-	-	
CD85k	ZM3.8	ASR	-	-	-	-	-	-	-	-	A46529	-	-	-	-	-	
CD86	HA5.2B7	ASR	-	-	-	-	IM2729U	-	-	B30647	B30648	-	-	-	-	B30646	
CD90	F15-42-1-5	ASR	-	-	IM1839U	-	IM1840U	-	-	-	-	-	-	-	-	-	
	Thy1/310	ASR	B30649	-	-	-	-	-	IM3703	-	-	-	-	-	-	B36121	

 New

AF 647 : Alexa Fluor* 647

AF 700 : Alexa Fluor* 700

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

ASR Single Color

			Laser		Blue Laser (488 nm)							Red Laser (640 nm)				
Emission Maximum			455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
Description	Clone	Status	PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD94	HP-3B1	ASR	B90465	-	-	-	-	-	-	-	-	B09980	-	-	-	-
CD99	HCD99	ASR	-	-	-	-	-	B76291	-	-	-	B76270	-	-	-	-
CD103	ZG5	ASR	-	-	IM1856U	-	-	-	-	-	-	B06204	-	-	-	-
CD105	TEA3/17.1.1	ASR	-	-	-	-	B76299	-	-	-	B43293	-	-	-	-	-
CD107a	H4A3	ASR	B13978	-	-	-	-	-	-	-	-	B12934	-	-	-	-
CD117	104D2D1	ASR	-	-	-	-	IM2732U	B38307	IM2733U	A66333	IM3698	IM3638	-	-	-	A86051
	95C3	ASR	-	-	-	-	IM1360U	-	IM2657U	-	-	-	-	-	-	-
CD122	CF1	ASR	-	-	-	-	-	-	-	-	A53365	-	-	-	-	-
CD123	SSDCLY107D2	ASR	-	-	-	-	A32535	B20027	-	B20022	B13647	B06376	-	-	B24028	-
	9F5	ASR	-	-	-	-	B14808	-	-	-	-	-	-	-	-	-
CD126	M91	ASR	-	-	-	-	IM1979	-	-	-	-	-	-	-	-	-
CD127	R34.34	ASR	-	-	B25366	-	IM1980U	-	A64617	-	A64618	B42026	-	-	A71116	B12700
CD133	W6B3C1	ASR	-	-	-	-	-	-	-	-	-	C15190	-	-	-	-
CD134	Ber-ACT35	ASR	-	-	-	-	-	-	-	-	-	B30629	-	-	-	-
CD135	SF1.340	ASR	-	-	-	-	IM2234U	-	-	-	-	-	-	-	-	-
CD137	4B4-1	ASR	-	-	-	-	-	B76262	-	-	-	-	-	-	-	-
CD138	B-A38	ASR	B37788	B42019	-	-	A40316	-	A40317	A98835	-	A87787	-	-	-	-
CD141	M80	ASR	-	-	B43289	-	-	-	-	-	-	-	-	-	-	-
CD144	TEA 1/31	ASR	-	-	-	-	A07481	-	-	-	-	-	-	-	-	-
CD146	TEA 1/34	ASR	-	-	-	-	A07483	-	A22364	-	-	-	-	-	-	-
CD154	TRAP-1	ASR	-	-	-	-	IM2216U	-	-	-	-	-	-	-	-	-
CD157	SY/11B5	ASR	-	-	-	-	B68160	-	-	-	-	-	-	-	-	-
CD158a,h	EB6B	ASR	-	-	-	-	A09778	-	-	A66898	A66899	A22332	-	-	-	-
CD158b1/b2,j	GL183	ASR	-	-	-	-	IM2278U	-	-	A66900	A66901	A22333	-	-	-	-
CD158e1/e2	Z27.3.7	ASR	-	-	-	-	IM3292	-	-	-	-	A60795	-	-	-	-
CD158f	UP-R1	ASR	-	-	-	-	B19715	-	-	-	-	B19716	-	-	-	-
CD158i	FES172	ASR	-	-	-	-	IM3337	-	-	-	-	A60796	-	-	-	-
CD159a	Z199	ASR	-	-	-	-	IM3291U	-	-	-	B10246	A60797	-	-	-	-
CD160	BY55	ASR	-	-	-	-	IM3657	-	-	-	-	-	-	-	-	-
CD161	191B8	ASR	-	-	-	-	IM3450	-	-	-	B30631	-	-	-	-	B30630
CD166	3A6	ASR	-	-	-	-	A22361	-	-	-	-	-	-	-	-	-
CD180	MHR73-11	ASR	-	-	-	-	B68148	-	-	-	-	-	-	-	-	-
CD183	G025H7	ASR	-	-	-	B68144	-	-	-	-	-	-	-	-	-	-
CD184	12G5	ASR	-	-	-	-	A07409	-	-	-	-	-	-	-	-	-
CD196	B-R35	ASR	-	-	-	-	-	-	-	-	B68132	-	-	-	-	-
CD197	G043H7	ASR	-	-	-	-	B30632	-	-	-	B46025	-	-	-	-	-
CD200	OX-104	ASR	-	-	-	-	B68124	-	-	-	B43299	-	-	-	-	B43301
CD203c	97A6	ASR	-	-	-	-	IM3575	-	-	-	A66906	-	-	-	-	-
CD206	3.29B1.10	ASR	B36119	-	-	-	IM2741	-	-	-	B36120	-	-	-	-	-
CD207	DCGM4	ASR	-	-	-	-	IM3577	-	-	-	-	-	-	-	-	-
CD235a	11E4B-7-6	ASR	-	-	IM2212U	-	IM2211U	-	-	B20026	A71564	-	-	-	-	A89314

 New

AF 647 : Alexa Fluor* 647

AF 700 : Alexa Fluor* 700

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

ASR Single Color

Description	Clone	Laser Emission Maximum	Status	Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)					
				455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
				PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD243	UIC2	ASR	-	-	-	-	IM2370U	-	-	-	-	-	-	-	-	-	-
CD244	C1.7	ASR	-	-	-	-	IM1608	-	-	B21171	-	-	-	-	-	-	-
CD274	PD-L1	ASR	-	-	-	-	-	-	-	-	-	B36126	-	-	-	-	-
CD278	ISA-3	ASR	-	-	-	-	-	-	-	-	-	B36127	-	-	-	-	-
CD279	PD1.3	ASR	-	-	-	-	B30634	-	-	B36123	A78885	B30633	-	-	-	-	-
CD294	BMI6	ASR	-	-	B42020	-	-	-	-	-	-	-	-	-	-	-	-
CD300a	E59.126	ASR	-	-	-	-	A22328	-	-	-	-	-	-	-	-	-	-
CD305	NKTA255	ASR	-	-	-	-	B68156	-	-	-	-	-	-	-	-	-	-
CD309	KDR-1	ASR	-	-	-	-	A64615	-	-	-	A64616	-	-	-	-	-	-
CD314	ON72	ASR	-	-	-	-	A08934	-	-	-	-	A22329	-	-	-	-	-
CD335	BAB281	ASR	-	-	-	-	IM3711	-	A66902	-	B38703	-	-	-	-	-	-
CD336	Z231	ASR	-	-	-	-	IM3710	-	A66903	-	-	-	-	-	-	-	-
CD337	Z25	ASR	-	-	-	-	IM3709	-	A66904	-	-	-	-	-	-	-	-
CD370	8F9	ASR	-	-	-	-	-	-	-	-	-	B43291	-	-	-	-	-
AP02.7	2.7A6A3	ASR	-	-	-	-	IM2088U	-	-	-	-	-	-	-	-	-	-
Cyclin A2	11B2G3	ASR	-	-	A22327	-	B15092	-	-	-	-	-	-	-	-	-	-
Cytokeratin	J1B3	ASR	-	-	IM2356U	-	-	-	-	-	-	-	-	-	-	-	-
FMC7	FMC7	ASR	B08466	-	IM1364U	-	-	-	-	-	-	-	-	-	-	-	-
FoxP3	259D	ASR	B90432	-	-	-	B46031	-	-	-	B46032	-	B30650	-	-	-	-
Granzyme B	GB11	ASR	-	-	B46038	-	-	-	-	-	-	-	-	-	-	-	-
HLA-ABC	B9.12.1	ASR	-	-	IM1838U	-	-	-	-	-	-	-	-	-	-	-	-
HLA-DR	Immu-357	ASR	A74781	B00070	IM1638U	-	IM1639U	IM3636	IM2659U	B20024	A40579	IM3635	-	-	-	-	B42021
	B8.12.2	ASR	-	-	IM0463U	-	IM0464U	-	-	-	-	-	-	-	-	-	-
HLA-DR, DP, DQ	9-49	ASR	-	-	6603424	-	6604366	-	-	-	-	-	-	-	-	-	-
IFN-γ	45,15	ASR	B46034	-	IM2716U	-	IM2717U	-	-	-	-	-	-	-	-	-	-
IL-2	N7.48A	ASR	-	-	-	-	IM2718U	-	-	-	-	-	-	-	-	-	-
IL-4	4D9	ASR	-	-	-	-	IM2719U	-	-	-	-	-	-	-	-	-	-
	MP4-25D2	ASR	-	-	-	-	-	-	-	-	B90441	-	-	-	-	-	-
Kappa	Rabbit polyclonal	ASR	-	-	A64828	-	-	-	-	-	-	-	-	-	-	-	-
	Goat polyclonal	ASR	-	-	C15623	-	-	-	-	-	-	B90420	-	-	-	-	-
Ki-67	Ki-67	ASR	-	-	-	B68180	-	-	-	-	-	-	-	-	-	-	-
Lambda	Rabbit polyclonal	ASR	-	-	-	-	A64827	-	-	-	-	-	-	-	-	-	-
	Goat polyclonal	ASR	-	-	-	-	C15189	-	-	-	B90416	-	-	-	-	-	-
Lactoferrin	CLB13.17	ASR	-	-	-	-	IM2742	-	-	-	-	-	-	-	-	-	-
Myeloperoxidase	CLB-MPO-1	ASR	-	-	IM1874U	-	IM3455U	-	-	B23132	-	-	-	-	-	-	-
NG2	7.1	ASR	-	-	-	-	IM3454U	-	-	-	-	-	-	-	-	-	-
NKp80	MA152	ASR	-	-	-	-	A52635	-	-	-	-	-	-	-	-	-	-
Perforin	dG9	ASR	B46030	-	-	-	-	-	-	-	-	-	-	-	-	-	-
ROR1	2A2	ASR	-	-	-	-	B68136	-	-	-	-	-	-	-	-	-	-
TCR PAN α/β	IP26A	ASR	-	-	-	-	A39499	-	A39500	-	-	B13981	-	-	-	-	-
TCR PAN γ/δ	IMMU 510	ASR	-	-	IM1571U	-	IM1418U	-	IM2662U	A99021	B10247	-	-	-	-	-	-

 New

AF 647 : Alexa Fluor* 647

AF 700 : Alexa Fluor* 700

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

ASR Single Color

		Laser	Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)					
		Emission Maximum	455 nm	528 nm	520 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm	780 nm
Description	Clone	Status	PB	KrO	FITC	AF 488	PE / RD1	ECD	PC5	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
TCR Vα24	C15	ASR	-	-	IM1589	-	IM2283	-	-	-	A66907	-	-	-	-	-
TCR Vβ1	BL37.2	ASR	-	-	IM2406	-	IM2355	-	-	-	-	-	-	-	-	-
TCR Vβ2	MPB2D5	ASR	-	-	IM2407	-	IM2213	-	-	-	-	-	-	-	-	-
TCR Vβ3	CH92	ASR	-	-	IM2372	-	-	-	-	-	-	-	-	-	-	-
TCR Vβ4	WJF24	ASR	-	-	B07084	-	IM3602	-	-	-	-	-	-	-	-	-
TCR Vβ5.1	IMMU 157	ASR	-	-	IM1552	-	IM2285	-	-	-	-	-	-	-	-	-
TCR Vβ5.2	36213	ASR	-	-	IM1482	-	IM2286	-	-	-	-	-	-	-	-	-
TCR Vβ5.3	3D11	ASR	-	-	-	-	IM2002	-	-	-	-	-	-	-	-	-
TCR Vβ7.1	ZOE	ASR	-	-	IM2408	-	IM2287	-	-	-	-	-	-	-	-	-
TCR Vβ7.2	ZIZOU4	ASR	-	-	B06666	-	IM3604	-	-	-	-	-	-	-	-	-
TCR Vβ8	56C5.2	ASR	-	-	IM1233	-	IM2289	-	-	-	-	-	-	-	-	-
TCR Vβ9	FIN9	ASR	-	-	-	-	IM2003	-	-	-	-	-	-	-	-	-
TCR Vβ11	C21	ASR	-	-	IM1586	-	IM2290	-	-	-	-	A66905	-	-	-	-
TCR Vβ12	VER2.32.1	ASR	-	-	IM1587	-	IM2291	-	-	-	-	-	-	-	-	-
TCR Vβ13.1	IMMU 222	ASR	-	-	IM1554	-	IM2292	-	-	-	-	-	-	-	-	-
TCR Vβ13.2	H132	ASR	-	-	-	-	IM3603	-	-	-	-	-	-	-	-	-
TCR Vβ13.6	JU74.3	ASR	-	-	IM1330	-	B07225	-	-	-	-	-	-	-	-	-
TCR Vβ14	CAS1.1.3	ASR	-	-	IM1558	-	IM2047	-	-	-	-	-	-	-	-	-
TCR Vβ16	TAMAYA1.2	ASR	-	-	IM1560	-	-	-	-	-	-	-	-	-	-	-
TCR Vβ17	E17.5F3.15.13	ASR	-	-	IM1234	-	IM2048	-	-	-	-	-	-	-	-	-
TCR Vβ18	BA62.6	ASR	-	-	-	-	IM2049	-	-	-	-	-	-	-	-	-
TCR Vβ20	ELL1.4	ASR	-	-	IM1562	-	IM2295	-	-	-	-	-	-	-	-	-
TCR Vβ21.3	IG125	ASR	-	-	IM1483	-	-	-	-	-	-	-	-	-	-	-
TCR Vβ22	IMMU 546	ASR	-	-	IM1484	-	IM2051	-	-	-	-	-	-	-	-	-
TCR Vβ23	AF23	ASR	-	-	-	-	IM2004	-	-	-	-	-	-	-	-	-
TCR Vδ1	R9.12	ASR	-	-	-	-	-	-	-	-	B49309	-	-	-	-	-
TCR Vδ2	IMMU 389	ASR	B49310	-	IM1464	-	-	-	-	-	-	-	-	-	-	-
TCR Vγ9	IMMU 360	ASR	-	-	IM1463	-	-	-	A63663	-	-	-	-	-	-	-
TCR ζ	2H2D9 (TIA-2)	ASR	-	-	-	-	IM3169	-	-	-	-	-	-	-	-	-
TdT	HTI+HT4+HT8+HT9	ASR	-	-	IM3524U	-	-	-	-	-	-	-	-	-	-	-
Thrombin receptor	SPAN 12	ASR	-	-	-	-	IM2583	-	-	-	-	-	-	-	-	-
	WEDE15	ASR	-	-	-	-	IM2584	-	-	-	-	-	-	-	-	-
TIA-1	2G9A10F5	ASR	-	-	-	-	IM3293	-	-	-	-	-	-	-	-	-
TNFα	IPM2 (188)	ASR	-	-	-	-	IM3279U	-	-	-	-	-	-	B76295	-	-
ZAP-70	SBZAP	ASR	-	-	-	-	B10248	-	-	-	-	-	-	-	-	-
ISOTYPIC CONTROLS																
IgG1 (mouse)	679.1Mc7	ASR	A74766	A96415	IM0639U	-	IM0670U	IM2714U	IM2663U	A62834	-	IM2475U	-	-	A71118	A71120
IgG2a (mouse)	7T4-1F5	ASR	-	-	A12690 B36624	-	A09141 B36633	A09144	A09147	-	A12692	A12694	-	-	-	-
IgG2b (mouse)	MPC-11	ASR	-	-	B36623	-	6603852	6604700	-	-	-	-	-	-	-	-
IgM (mouse)	GC323	ASR	-	-	IM1269U	-	-	-	-	-	-	-	-	-	-	-
	R4A3-22-12	ASR	-	-	B36627	-	-	-	-	-	-	-	-	-	-	-

 New

AF 647 : Alexa Fluor® 647

AF 700 : Alexa Fluor® 700

APC-A700 : APC-Alexa Fluor® 700

APC-A750 : APC-Alexa Fluor® 750

PB : Pacific Blue®

KrO : Krome Orange

IVD Single Color

Description	Clone	Status	Laser		Blue Laser (488 nm)						Red Laser (640 nm)				
			Emission Maximum	455 nm	528 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	668 nm	719 nm	719 nm
			PB	KrO	FITC	PE / RD1	ECD	PCS	PC5.5	PC7	APC	AF 647	AF 700	APC-A700	APC-A750
CD2	SFC13Pt2H9 (T11)	IVD & CE	-	-	6602389	6602868	-	-	-	-	-	-	-	-	-
CD3	UCHT1	IVD & CE	-	-	6604623 6604625	6604621 6604627	-	-	-	-	-	-	-	-	-
CD4	SFC112T4D11 (T4)	IVD & CE	-	-	6603862 6602393	6603850 6602864	-	-	-	-	-	-	-	-	-
CD5	SFC124T6G12 (T1)	IVD & CE	-	-	-	6603020	-	-	-	-	-	-	-	-	-
CD8	SFC121Thy2D3 (T8)	IVD & CE	-	-	6603861 6602385	6603848 6603274	-	-	-	-	-	-	-	-	-
CD14	322A-1	IVD & CE	-	-	6603511	6603262	-	-	-	-	-	-	-	-	-
	T16	IVD	-	-	6602507	6603266	-	-	-	-	-	-	-	-	-
CD19	89B	IVD & CE	-	-	6603012	6603024	-	-	-	-	-	-	-	-	-
CD20	H299	IVD & CE	-	-	6602381	6603446	-	-	-	-	-	-	-	-	-
ISOTYPIC CONTROLS															
IgG1 (mouse)	ZT8-2F5	IVD & CE	-	-	6602928 6603864	6602884	-	-	-	-	-	-	-	-	-

ASR : Analyte Specific Reagents. Analytical and performance characteristics are not established.

IVD: In Vitro Diagnostic Product.

IVD & CE: In Vitro Diagnostic Products & CE-Marked Products.

AF 647 : Alexa Fluor® 647

AF 700 : Alexa Fluor® 700

APC-A700 : APC-Alexa Fluor® 700

APC-A750 : APC-Alexa Fluor® 750

PB : Pacific Blue*

KrO : Krome Orange

 New

* Alexa Fluor and Pacific Blue are registered trademarks of Molecular Probes, Inc.

Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

FLOW-4018SBO9.18

© 2018 Beckman Coulter, Inc.

PRINTED IN U.S.A.

