



CE-IVD SINGLE COLOR ANTIBODIES

OUS Version



CE-IVD Single Color

Description	Clone	Status	Laser		Emission Maximum						Laser		
			Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)		
			455 nm	528 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	719 nm	780 nm
			PB	KrO	FITC	PE / RD1	ECD	PC5	PC5.5	PC7	APC	APC-A700	APC-A750
ANTIBODIES TO CD ANTIGENS													
CD1a	BL6	CE	-	-	-	A07742	-	-	-	-	-	-	-
CD2	39C1.5	CE	-	-	A07743	A07744	-	A07745	-	A21689	B49210	-	-
	SFC13Pt2H9	IVD & CE	-	-	6602389 6603863	6602868 6603849	-	-	-	-	-	-	-
CD3	UCHT1	CE	B49204	-	A07746	A07747	A07748	A07749	B49203	737657	IM2467	-	A94680
		IVD & CE	-	-	6604623 6604625	6604621 6604627	-	-	-	-	-	-	-
CD4	13B8.2	CE	B49197	-	A07750	A07751	-	A07752	-	-	IM2468	-	A94682
	SFC112T4D11	IVD & CE	-	-	6602393 6603862	6602864 6603850	-	-	-	-	-	-	-
		CE	-	-	-	-	-	-	-	737660	-	-	-
CD5	BL1a	CE	-	-	A08932	A07753	-	A07754	B49191	A21690	-	-	B92413
	SFC124T6G12	IVD & CE	-	-	-	6603020	-	-	-	-	-	-	-
CD7	8H8.1	CE	-	-	A07755	-	-	IM3613	-	B49183	-	B36290	-
CD8	B9.11	CE	B49182	-	A07756	A07757	-	A07758	-	-	IM2469	B49181	A94683
	SFC121Thy2D3	IVD & CE	-	-	6602385 6603861	6603274 6603848	-	-	-	-	-	-	-
		CE	-	-	-	-	737659	-	-	737661	-	-	-
CD10	ALB1	CE	-	-	A07759	A07760	-	A07761	-	B96750	B92400	B49223	-
CD11b	Bear1	CE	B36296	-	IM0530	-	-	-	-	-	-	-	B36295
CD11c	BU15	CE	-	-	-	IM1760	-	-	-	B96763	-	-	-
CD13	SJ1D1	CE	-	-	-	A07762	-	-	-	-	-	-	-
	Immu103.44	CE	-	-	-	-	B36286	A07763	B49196	-	-	-	-
CD14	RM052	CE	-	-	B36297	A07764	B92391	A07765	-	-	IM2580	-	B92421
	322A-1	IVD & CE	-	-	6603511 6604110	6603262	-	-	-	-	-	-	-
	116	IVD & CE	-	-	6604114	6604500	-	-	-	-	-	-	-
IVD		-	-	6602507	6603266	-	-	-	-	-	-	-	
CD15	80H5	CE	B49218	-	B36298	-	-	B49217	-	-	-	-	-
CD16	3G8	CE	B36292	-	B49215	A07766	B49216	A07767	-	-	-	-	B49184
CD19	J3-119	CE	B49213	-	A07768	A07769	A07770	A07771	B49211	IM3628	IM2470	B49212	A94681
	89B	IVD & CE	-	-	6603012 6603859	6603024 6603846	-	-	-	-	-	-	-
CD20	B9E9	CE	B49208	-	A07772	IM1451	B92433	A07773	-	IM3629	A21693	-	B49209
	H299	IVD & CE	-	-	6602381 6604106	6603446 6603858	-	-	-	-	-	-	-
	FMC7	CE	B96782	-	A07791	-	-	-	-	-	-	-	-
CD22	SJ10.1H11	CE	-	-	-	IM1835	-	-	-	-	B96777	B36293	-
CD23	9P25	CE	-	-	IM0529	B49207	-	-	-	-	-	-	-
CD24	ALB9	CE	-	-	-	B92425	-	-	-	-	-	-	-

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

CE-IVD Single Color

Description	Clone	Status	Laser		Emission Maximum						Laser		
			Violet Laser (405 nm)		Blue Laser (488 nm)						Red Laser (640 nm)		
			455 nm	528 nm	520 nm	575 nm	620 nm	670 nm	694 nm	767 nm	660 nm	719 nm	780 nm
			PB	KrO	FITC	PE / RD1	ECD	PC5	PC5.5	PC7	APC	APC-A700	APC-A750
CD25	B1.49.9	CE	-	-	-	A07774	-	IM2646	B92458	-	-	B92454	-
CD27	1A4CD27	CE	-	-	-	B96790	-	-	-	B49205	-	-	-
CD33	D3HL60.251	CE	-	-	-	A07775	-	IM2647	B36289	B92408	IM2471	-	-
CD34	581	CE	-	-	IM1870	A07776	B49202	A07777	-	A21691	IM2472	B92417	B92463
CD36	FA6.152	CE	-	-	B49201	-	-	-	-	-	-	-	-
CD38	LS198-4-3	CE	B92396	-	-	A07779	-	A07780	B49199	B49198	-	-	B49200
	T16	CE	-	-	A07778	-	-	-	-	-	-	-	-
CD41	P2	CE	-	-	-	A07781	-	-	-	-	-	-	-
CD43	DFT1	CE	-	-	-	-	-	-	-	-	-	-	B49195
CD45	J33	CE	A74763	B36294	A07782	A07783	A07784	A07785	A62835	IM3548	IM2473	A79390	A79392
CD45RA	ALB11	CE	-	-	A07786	-	-	-	-	-	-	-	-
	ZH4LDH11LDB9	CE	-	-	-	-	B49193	-	-	-	-	-	B49194
CD45RO	UCHL1	CE	-	-	-	A07787	B49192	-	-	-	-	-	-
CD55	JS11KSC2.3	CE	-	-	-	B49190	-	-	-	-	-	-	-
CD56	N901	CE	-	-	-	A07788	B49214	A07789	B49189	A21692	IM2474	B92446	-
CD57	NC1	CE	-	-	B49188	-	-	-	-	-	-	-	-
CD59	P282E	CE	-	-	B49187	-	-	-	-	-	-	-	-
CD61	SZ21	CE	-	-	IM1758	-	-	-	-	-	-	-	-
CD62P	CLB-Thromb/6	CE	-	-	A07790	-	-	-	-	-	-	-	-
CD63	CLB-Grand/12	CE	-	-	B92467	-	-	-	-	-	-	-	-
CD64	22	CE	-	-	B49185	-	-	B49186	-	-	-	-	B96769
CD65	88H7	CE	-	-	B36299	-	-	-	-	-	-	-	-
CD69	TP1.55.3	CE	-	-	-	-	-	IM2656	-	-	-	-	-
CD71	YDJ1.2.2	CE	-	-	IM0483	-	-	-	-	-	-	-	-
CD79a	HM47	CE	-	-	-	IM2221	-	-	-	-	B36287	-	-
CD79b	CB3-1	CE	-	-	-	IM1612	-	-	-	-	B96773	-	-
CD103	2G5	CE	-	-	B49222	-	-	-	-	-	-	-	-
CD105	TEA3/17.1.1	CE	-	-	-	B92442	-	-	-	-	-	-	-
CD117	104D2D1	CE	-	-	-	IM2732	-	IM2733	B96754	B49221	B36300	-	B92450
CD127	R34.34	CE	-	-	-	B49220	-	-	-	-	-	-	-
CD138	B-A38	CE	-	-	-	A54190	-	A54191	B96786	-	B49219	-	-
CD200	OX-104	CE	-	-	-	-	-	-	-	B92472	-	-	-
CD203c	97A6	CE	-	-	-	B92404	-	-	-	-	-	-	-
CD235a	11E4B-7-6	CE	-	-	B49206	A07792	-	-	-	-	-	-	-

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

CE-IVD Single Color

		Laser	Violet Laser (405 nm)		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	719 nm	780 nm
Description	Clone	Status	PB	KrO	FITC	PE / RD1	ECD	PC5	PC5.5	PC7	APC	APC-A700	APC-A750
ANTIBODIES TO NON CD ANTIGENS													
FMC7	FMC7	CE	B96782	-	A07791	-	-	-	-	-	-	-	-
HLA-DR	Immu-357	CE	B36291	-	B96758	IM1639	B92438	A07793	-	B49180	-	-	-
κ chain	Polyclonal	CE	-	-	B49178	-	-	-	-	-	-	-	-
λ chain	Polyclonal	CE	-	-	-	B49174	-	-	-	-	-	-	-
Myeloperoxidase	CLB-MPO-1	CE	-	-	IM1874	B36288	-	-	-	-	-	-	-
NG2	7.1	CE	-	-	-	B92429	-	-	-	-	-	-	-
TCR PAN α/β	IP26A	CE	-	-	-	B49177	-	-	-	-	-	-	-
TCR PAN γ/δ	IMMU510	CE	-	-	B49175	B49176	-	IM2662	-	-	-	-	-
TdT	HT1 + HT4 + HT8 + HT9	CE	-	-	IM3524	-	-	-	-	-	-	-	-
ISOTYPIC CONTROLS													
IgG1 (Isotype)	679.1Mc7	CE	A74764	-	A07795	A07796	A07797	A07798	A62833	737662	IM2475	A79391	A79393
	2T8-2F5	IVD & CE	-	-	6602928 6603864	6602884	-	-	-	-	-	-	-
IgG2a (Isotype)	7T4-1F5	CE	-	-	A12689	A09142	-	A09148	-	-	A12693	-	-
		IVD & CE	-	-	6603855	6604119	-	-	-	-	-	-	-
IgG2b (Isotype)	MPC-11	IVD & CE	-	-	6603853	-	-	-	-	-	-	-	-

CE: In vitro diagnostic products compliant with EU directive 98/79/EC.

IVD: In Vitro Diagnostic products.

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

CE-marked reagents are sold in European Union countries and in countries where they are compliant with local regulation (not in United States).

IVD products are sold in countries where they comply with applicable regulation.

APC-A700 : APC-Alexa Fluor* 700

APC-A750 : APC-Alexa Fluor* 750

PB : Pacific Blue *

KrO : Krome Orange

* 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-4017SB09.18

© 2018 Beckman Coulter, Inc.

