



Miltenyi Biotec

REAscreen™ MAX Kit, human, FFPE, version 01

Panel composition – 89 antibody conjugates

130-127-847

REAscreen MAX Antibody Plate

	1	2	3	4	5	6	7	8	9	10	11	12
A	CD158i (KIR2DS4)	IRF-7 pS477/pS479	SSRP1	CD45RA	ZAP70	CD181 (CXCR1)	TTF-1	CD138	Sox2	CD104 (Integrin β4)	CD295 (LEPR)	CD162
B	CD134 (OX40)	CD240DCE	IRF-7	CD15	CD43	AN2	Galectin-3	CD66c	TOM22	CD147	CD66acde	
C	Syk	Slug	LRP-4	IgD	Cyto-keratin	Vimentin	CLA	CD279 (PD1)	Dectin-1	Galectin-9	CD234 (DARC)	
D	PAX-5 (BSAP)	FAK pS910	p53	FcεR1α	PCNA	CD195 (CCR5)	CD268	CD196 (CCR6)	CD88 (C5AR)	BMI-1	CD305 (LAIR-1)	
E	PKCa	BATF	CD334 (FGFR4)	CD45R (B220)	Bcl-2	CD226 (DNAM-1)	CD65	CD223	CD52	CD66abce	AKT Pan (PKB)	
F	CD182 (CXCR2)	CD20 Cyto-plasmic	CD45	CD326 (EpCAM)	Cyto-keratin 8	Podoplanin	MLC2v	HLA-DR, DP, DQ	CD66b	Hsp70	CD45RB	
G	CD202b (TIE-2)	Dnmt3b	Jak1	SUSD2	EZH2	CD235a (Glycophorin A)	α-Actinin (Sarcomeric)	CD79a	CD40	CD99	CD317 (BST2)	
H	CD239 (BCAM)	p120 Catenin pS879	CD44	CD57	CD123	Ki-67	CD233	CD45RO	β-Catenin	CD271 (LNGFR)	CD100	

Antigen	Clone	Fluorochrome	Antibody type	Plate position
AKT Pan (PKB)	REA676	PE	REAffinity™	E11
AN2	REA989	PE	REAffinity	B6
BATF	REA486	FITC	REAffinity	E2

Antigen	Clone	Fluorochrome	Antibody type	Plate position
Bcl-2	REA872	FITC	REAffinity	E5
BMI-1	REA438	FITC	REAffinity	D10
CD100	REA316	FITC	REAffinity	H11
CD104 (Integrin β 4)	REA236	FITC	REAffinity	A10
CD123	REA918	FITC	REAffinity	H5
CD134 (OX40)	ACT35	FITC	Hybridoma	B1
CD138	44F9	FITC	Hybridoma	A8
CD147	REA282	FITC	REAffinity	B10
CD15	VIMC6	FITC	Hybridoma	B4
CD158i (KIR2DS4)	JJC11.6	PE	Hybridoma	A1
CD162	REA319	FITC	REAffinity	A12
CD181 (CXCR1)	REA958	FITC	REAffinity	A6
CD182 (CXCR2)	REA208	PE	REAffinity	F1
CD195 (CCR5)	REA245	PE	REAffinity	D6
CD196 (CCR6)	REA190	PE	REAffinity	D8
CD20 Cytoplasmic	REA543	PE	REAffinity	F2
CD202b (TIE-2)	REA198	PE	REAffinity	G1
CD223	REA351	PE	REAffinity	E8
CD226 (DNAM-1)	REA1040	FITC	REAffinity	E6
CD233	REA368	FITC	REAffinity	H7
CD234 (DARC)	REA376	PE	REAffinity	C11
CD235a (Glycophorin A)	REA175	FITC	REAffinity	G6
CD239 (BCAM)	REA276	PE	REAffinity	H1
CD240DCE	REA327	PE	REAffinity	B2
CD268	REA1115	FITC	REAffinity	D7
CD271 (LNGFR)	REAL709	APC	REAlEase®	H10
CD279 (PD1)	REA1165	PE	REAffinity	C8
CD295 (LEPR)	REA361	PE	REAffinity	A11
CD305 (LAIR-1)	REA447	PE	REAffinity	D11
CD317 (BST2)	REA202	FITC	REAffinity	G11
CD326 (EpCAM)	REA764	FITC	REAffinity	F4
CD334 (FGFR4)	4FR6D3	FITC	Hybridoma	E3
CD40	HB14	PE	Hybridoma	G9
CD43	REA833	FITC	REAffinity	B5
CD44	REA690	FITC	REAffinity	H3
CD45	5B1	FITC	Hybridoma	F3
CD45R (B220)	REA755	FITC	REAffinity	E4
CD45RA	REA562	PE	REAffinity	A4
CD45RB	REA119	FITC	REAffinity	F11
CD45RO	REAL377	PE	REAlEase	H8
CD52	REA164	FITC	REAffinity	E9
CD57	REA769	PE	REAffinity	H4
CD65	REA1128	PE	REAffinity	E7
CD66abce	REA1230	PE	REAffinity	E10
CD66acde	REA428	PE	REAffinity	B11
CD66b	REA306	FITC	REAffinity	F9
CD66c	REA414	FITC	REAffinity	B8
CD79a	REA1168	PE	REAffinity	G8
CD88 (C5AR)	REA1213	PE	REAffinity	D9
CD99	REAL687	PE	REAlEase	G10
CLA	REA1101	PE	REAffinity	C7
Cytokeratin	REA831	FITC	REAffinity	C5

Antigen	Clone	Fluorochrome	Antibody type	Plate position
Cytokeratin 8	REA885	FITC	REAffinity	F5
Dectin-1	REA515	PE	REAffinity	C9
Dnmt3b	REA565	PE	REAffinity	G2
EZH2	REA907	FITC	REAffinity	G5
FAK pS910	REA407	PE	REAffinity	D2
FcεR1α	REA758	FITC	REAffinity	D4
Galectin-3	REA1072	FITC	REAffinity	B7
Galectin-9	REA435	FITC	REAffinity	C10
HLA-DR, DP, DQ	REA332	PE	REAffinity	F8
Hsp70	REA349	FITC	REAffinity	F10
IgD	REA740	PE	REAffinity	C4
IRF-7	REA521	FITC	REAffinity	B3
IRF-7 pS477/pS479	REA310	PE	REAffinity	A2
Jak1	REA700	PE	REAffinity	G3
Ki-67	REA183	FITC	REAffinity	H6
LRP-4	REA552	PE	REAffinity	C3
MLC2v	REA401	PE	REAffinity	F7
p120 Catenin pS879	REA548	PE	REAffinity	H2
p53	REA609	FITC	REAffinity	D3
PAX-5 (BSAP)	REA140	PE	REAffinity	D1
PCNA	REA858	FITC	REAffinity	D5
PKCα	REA156	PE	REAffinity	E1
Podoplanin	REA446	PE	REAffinity	F6
Slug	REA404	PE	REAffinity	C2
Sox2	REA320	PE	REAffinity	A9
SSRP1	REA555	FITC	REAffinity	A3
SUSD2	REA795	PE	REAffinity	G4
Syk	REA111	FITC	REAffinity	C1
TOM22	REA1185	PE	REAffinity	B9
TTF-1	REA1090	FITC	REAffinity	A7
Vimentin	REA409	FITC	REAffinity	C6
ZAP70	REA814	PE	REAffinity	A5
α-Actinin (Sarcomeric)	REA402	FITC	REAffinity	G7
β-Catenin	REA480	FITC	REAffinity	H9



Miltenyi Biotec B.V. & Co. KG | Phone +49 2204 8306-0 | Fax +49 2204 85197 | macsde@miltenyi.com | www.miltenyibiotec.com

Miltenyi Biotec provides products and services worldwide. Visit www.miltenyibiotec.com/local to find your nearest Miltenyi Biotec contact.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. MACS and the Miltenyi Biotec logo are registered trademarks or trademarks of Miltenyi Biotec and/or its affiliates in various countries worldwide. Copyright © 2021 Miltenyi Biotec and/or its affiliates. All rights reserved.