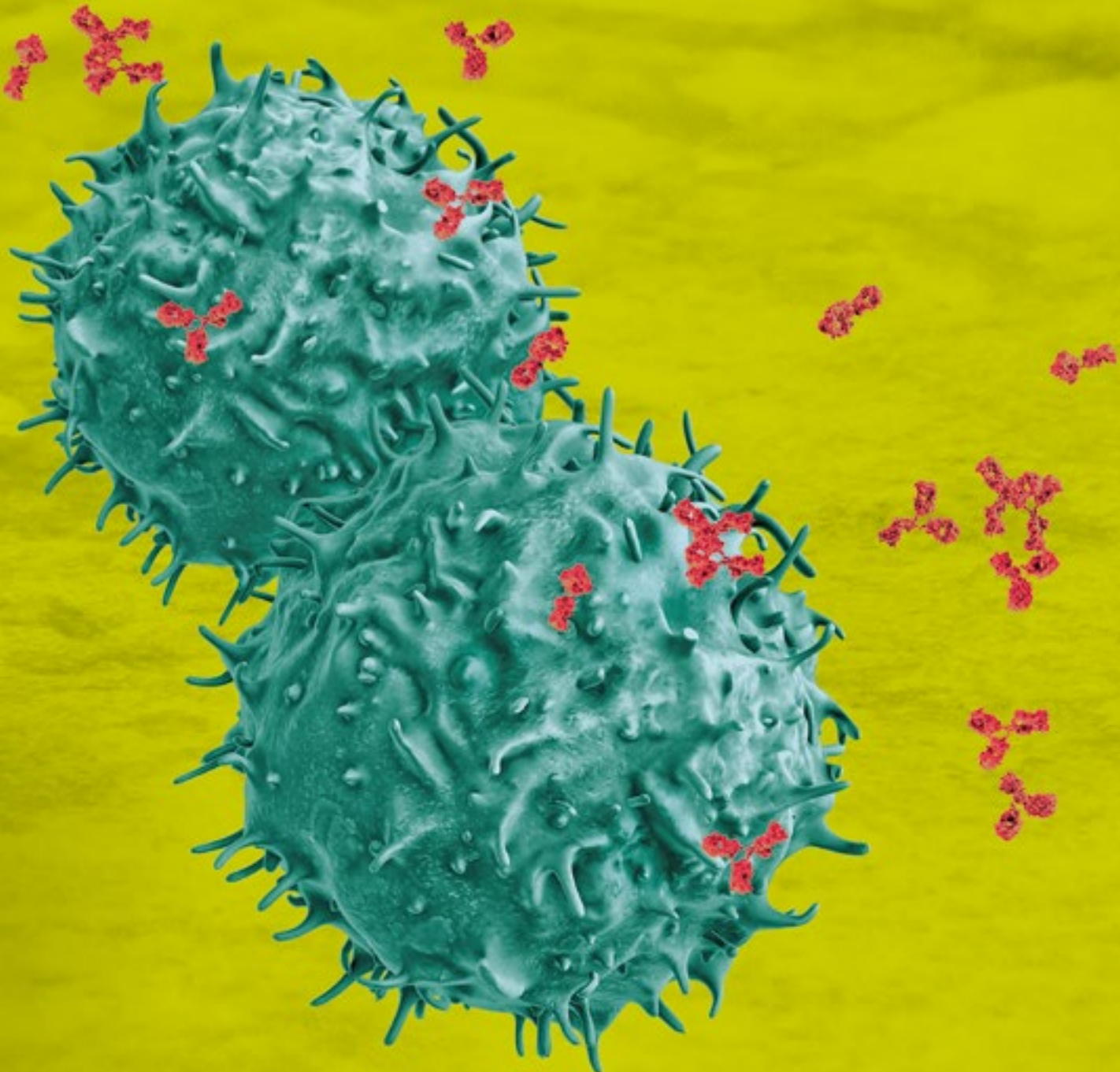




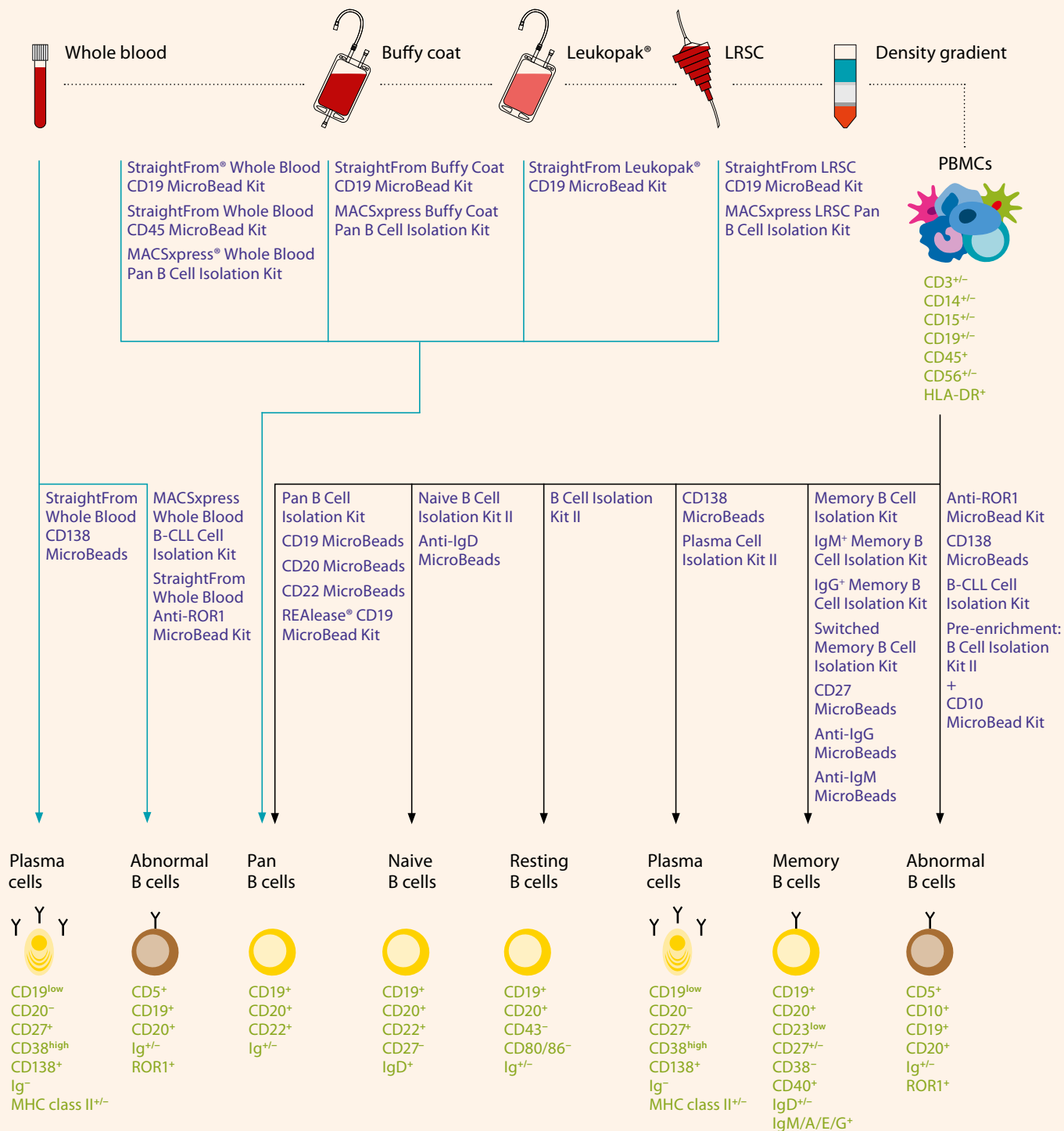
Miltenyi Biotec

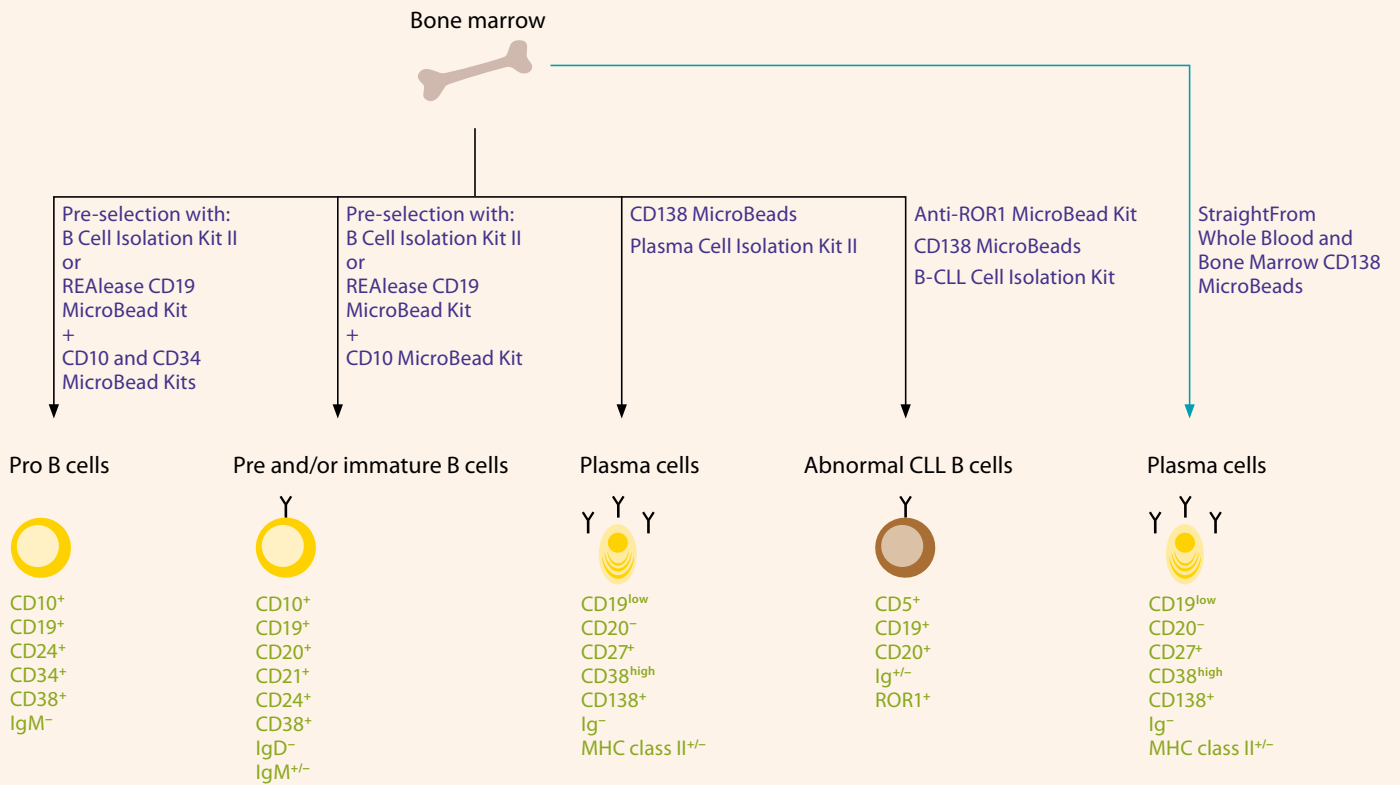
The complete portfolio for your B cell research

Product quick guide



Human B cells from whole blood and bone marrow





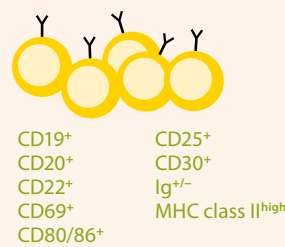
B cell cultivation and activation

B cells and B cell subsets



B Cell Expansion Kit, TLR-Ligands*
Cytokines, e.g., IL-4*, IL-21, TNF-α*
Human CD40-Ligand Multimer Kits

Expanded and activated B cells



Legend

- direct cell separation products
- separation
- differentiation/cultivation

- MACS[®] Cell Separation products (RUO)
- MACS Cell Culture and Stimulation products
- MACS Flow Cytometry and Cytokine detection products (RUO)

- CD:** cluster of differentiation
- CLL:** chronic lymphocytic leukemia
- HLA:** human leukocyte antigen
- Ig:** immunoglobulin
- LRSC:** leukocyte reduction system chamber
- MHC:** major histocompatibility complex
- PBMC:** peripheral blood mononuclear cells
- ROR:** receptor tyrosine kinase-like orphan receptor 1
- TNF:** tumor necrosis factor

- no expression
- + expression
- ^{low} low expression
- ^{int} intermediate expression
- ^{high} high expression
- +/- expression dependent on subset

*Available also as MACS GMP Product

Human B cell research

Product tables

Cell separation reagents

Products	Order no.
Anti-IgD MicroBeads	130-103-775
Anti-IgG MicroBeads	130-047-501
Anti-IgM MicroBeads	130-093-230
Anti-ROR1 MicroBead Kit	130-103-929
B Cell Isolation Kit II	130-091-151
B-CLL Cell Isolation Kit	130-103-466
CD10 MicroBead Kit	130-093-452
CD19 MicroBeads	130-050-301
CD20 MicroBeads	130-091-104
CD22 MicroBeads	130-046-401
CD27 MicroBeads	130-051-601
CD38 MicroBead Kit	130-092-263
CD43 MicroBeads	130-091-333
CD138 MicroBeads	130-051-301
IgG ⁺ Memory B Cell Isolation Kit	130-094-350
IgM ⁺ Memory B Cell Isolation Kit	130-093-619
Memory B Cell Isolation Kit	130-093-546
Naive B Cell Isolation Kit II	130-091-150
Pan B Cell Isolation Kit	130-101-638
Plasma Cell Isolation Kit II	130-093-628
Switched Memory B Cell Isolation Kit	130-093-617
REAl ease Technology	
REAl ease CD19 MicroBead Kit	130-117-034
StraightFrom Technology	
StraightFrom Buffy Coat CD19 MicroBead Kit	130-114-974
StraightFrom Leukopak [®] CD19 MicroBead Kit	130-117-021
StraightFrom LRSC CD19 MicroBead Kit	130-117-025
StraightFrom Whole Blood CD19 MicroBeads	130-090-880
StraightFrom Whole Blood CD138 MicroBeads	130-093-062
StraightFrom Whole Blood and Bone Marrow CD138 MicroBeads	130-105-961
StraightFrom Whole Blood Anti-ROR1 MicroBead Kit	130-105-751
MACSxpress Technology	
MACSxpress Buffy Coat Pan B Cell Isolation Kit	130-120-002
MACSxpress LRSC Pan B Cell Isolation Kit	130-119-998
MACSxpress Whole Blood B-CLL Cell Isolation Kit	130-104-445
MACSxpress Whole Blood Pan B Cell Isolation Kit	130-104-444

Cell culture reagents

Products	Research grade	Premium grade	MACS GMP
BAFF	●		
B Cell Expansion Kit		●	
CD40-Ligand Multimer Kit		●	
CXCL13	●		
Flt3-Ligand	●	●	●
IFN-γ1b	●	●	
IL1-beta	●	●	●
IL-2 improved sequence (IS)	●	●	●
IL-3	●	●	●
IL-4	●	●	●
IL-6	●	●	●
IL-7	●	●	●
IL-10	●		
IL-12		●	
IL-13	●	●	
IL-17	●		
IL-21	●	●	●
IL-35	●		
PepTivator [®] Peptide Pools	●	●	●
SCF	●	●	●
SDF-1α	●		
TGF-β1		●	●
TLR7-Ligands	●		
TLR9-Ligands	●		●
TNF-α	●	●	●
TSLP	●		

For more information about custom fillings and cocktails visit:

www.miltenyibiotec.com/cellculture

Flow cytometry antibodies

Antigen	Hybridoma Clones	REAffinity™ Clones	REALease Clones
CD1d	5I.1		
CD5	UCHT2	REA782	REAL382
CD10	97C5	REA877	REAL318
CD19	LT19	REA675	REAL106
CD20	LT20	REA780	
CD21	HB5	REA940	
CD22		REA340	
CD23	M-L23.4	REA1222	
CD24	32D12	REA832	
CD25	4E3, 3G10	REA570, REA945	REAL128
CD27	M-T271, LG.3A10	REA499	REAL173
CD32	2E1	REA997	
CD38	IB6	REA572, REA671	
CD43	DF-T1	REA833	
CD45	5B1	REA747	REAL258
CD45RA	T6D11	REA562, REA1047	
CD45RB		REA119	
CD45RO	UCLH1	REA611	REAL377
CD80	2D10	REA661	
CD86	FM95	REA968	
CD93	VIMD2	REA1111	
CD95 (FAS)	DX2	REA738	REAL186
CD138	44F9		
IL-10	JES3-9D7	REA842	
IgD	IgD26	REA740	
IgM	PJ2-22H3		
IgG	IS11-3B2.2.3		
IgG (H+L)	Polyclonal		
IgG1	IS11-12E4.23.20		
IgE	MB10-5C4	REA1049	
IgA	IS11-8E10	REA1014	
IgA2	IS11-21E11	REA995	
HLA	Many different clones to several HLA class antigens		

Flow cytometry cocktails and kits

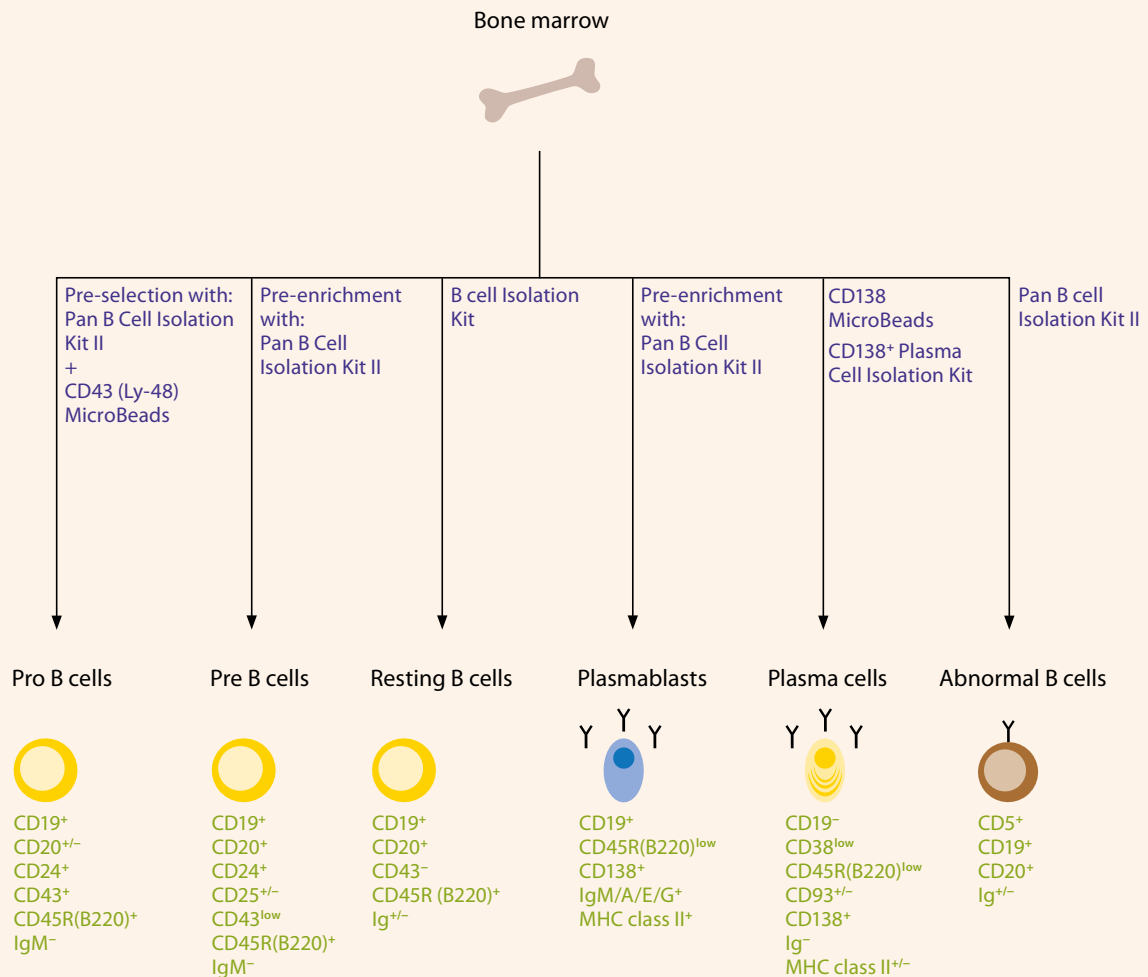
Products	Order no.
GM-CSF Secretion Assay – Detection Kit (PE)	130-105-755
IFN-γ Secretion Assay – Detection Kit (APC)	130-090-762
IFN-γ Secretion Assay – Detection Kit (FITC)	130-090-433
IFN-γ Secretion Assay – Detection Kit (PE)	130-054-202
IL-2 Secretion Assay – Detection Kit (APC)	130-090-763
IL-2 Secretion Assay – Detection Kit (PE)	130-090-487
IL-4 Secretion Assay – Detection Kit (PE)	130-054-102
IL-5 Secretion Assay – Detection Kit (APC)	130-091-624
IL-5 Secretion Assay – Detection Kit (PE)	130-091-623
IL-10 Secretion Assay – Detection Kit (APC)	130-090-761
IL-10 Secretion Assay – Detection Kit (PE)	130-090-434
IL-12 Secretion Assay – Cell Enrichment and Detection Kit (PE)	130-092-122
RCI Anti-IL-2-PE	130-099-450
RCI Anti-IFN-γ-APC	130-097-596
RCI Anti-IFN-γ-PE	130-097-600
TNF-α Secretion Assay – Detection Kit (APC)	130-091-267
TNF-α Secretion Assay – Detection Kit (PE)	130-091-268

MACSPlex Custom Cytokine Assays

Products	Order no.
MACSPlex Cytokine Basic Kit, human and mouse	130-109-701
MACSPlex Cytokine 12 Kit	130-099-169
MACSPlex GM-CSF Reagents Kit	130-109-684
MACSPlex IFN-α Reagents Kit	130-109-683
MACSPlex IFN-γ Reagents Kit	130-109-695
MACSPlex IL-2 Reagents Kit	130-109-704
MACSPlex IL-4 Reagents Kit	130-109-689
MACSPlex IL-5 Reagents Kit	130-109-705
MACSPlex IL-6 Reagents Kit	130-109-691
MACSPlex IL-9 Reagents Kit	130-109-693
MACSPlex IL-10 Reagents Kit	130-109-699
MACSPlex IL-12 Reagents Kit	130-109-697
MACSPlex IL-17A Reagents Kit	130-109-688
MACSPlex TNF-α Reagents Kit	130-109-694

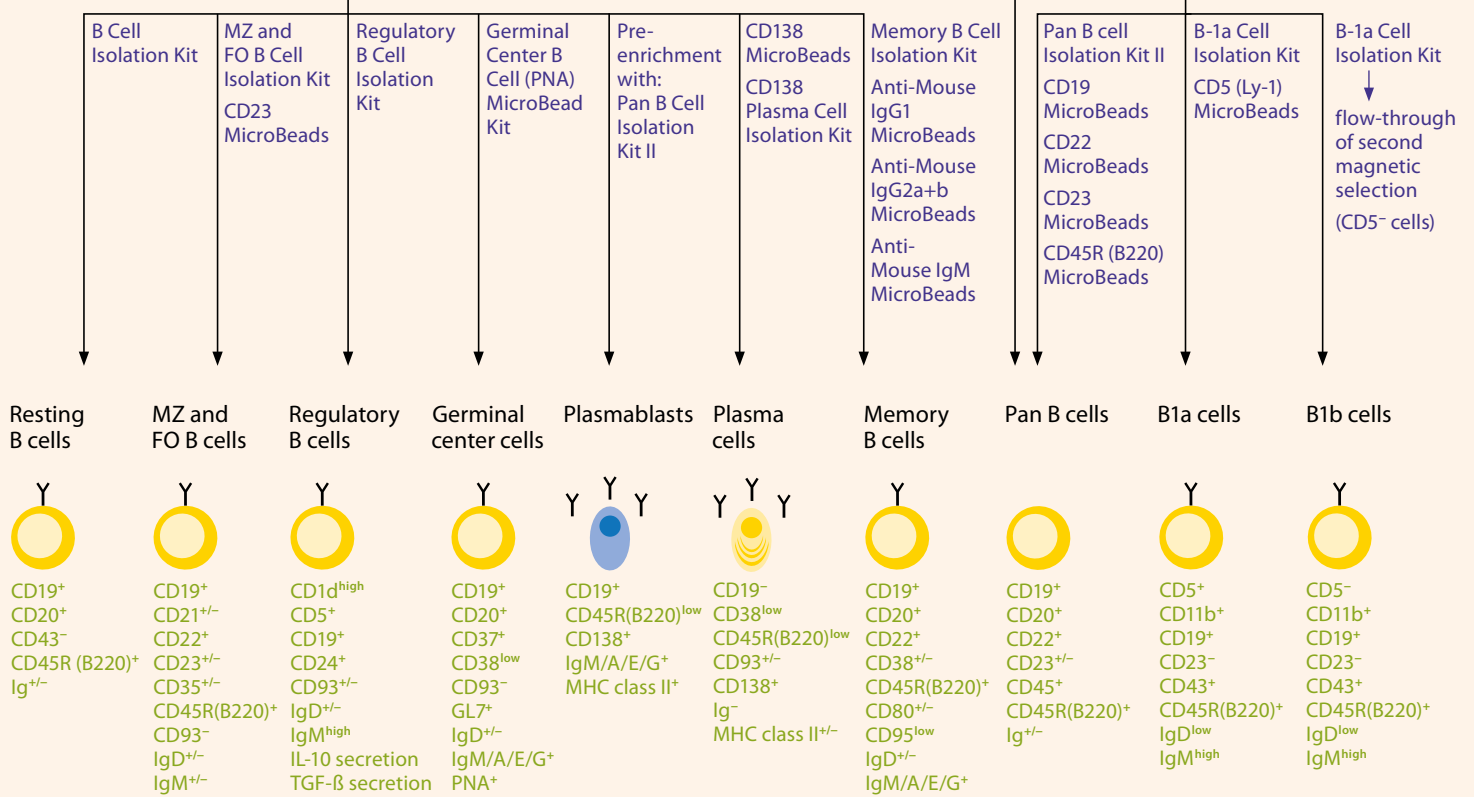
For more information about antibodies and fluorescent conjugates for flow cytometry visit:
www.miltenyibiotec.com/flowcytometry

Mouse B cells from bone marrow, spleen, and lymphoid tissue



Spleen/lymphoid tissue

Body cavities



Legend

→ separation

■ MACS Cell Separation products (RUO)

■ MACS Flow Cytometry and Cytokine Detection products (RUO)

CD: cluster of differentiation

FO: follicular

Ig: immunoglobulin

MHC: major histocompatibility complex

MZ: marginal zone

PNA: peanut agglutinin

TGF: transforming growth factor

- no expression

^{low} low expression

^{int} intermediate expression

+ expression

^{high} high expression

^{+/-} expression dependent on subset

Mouse B cell research

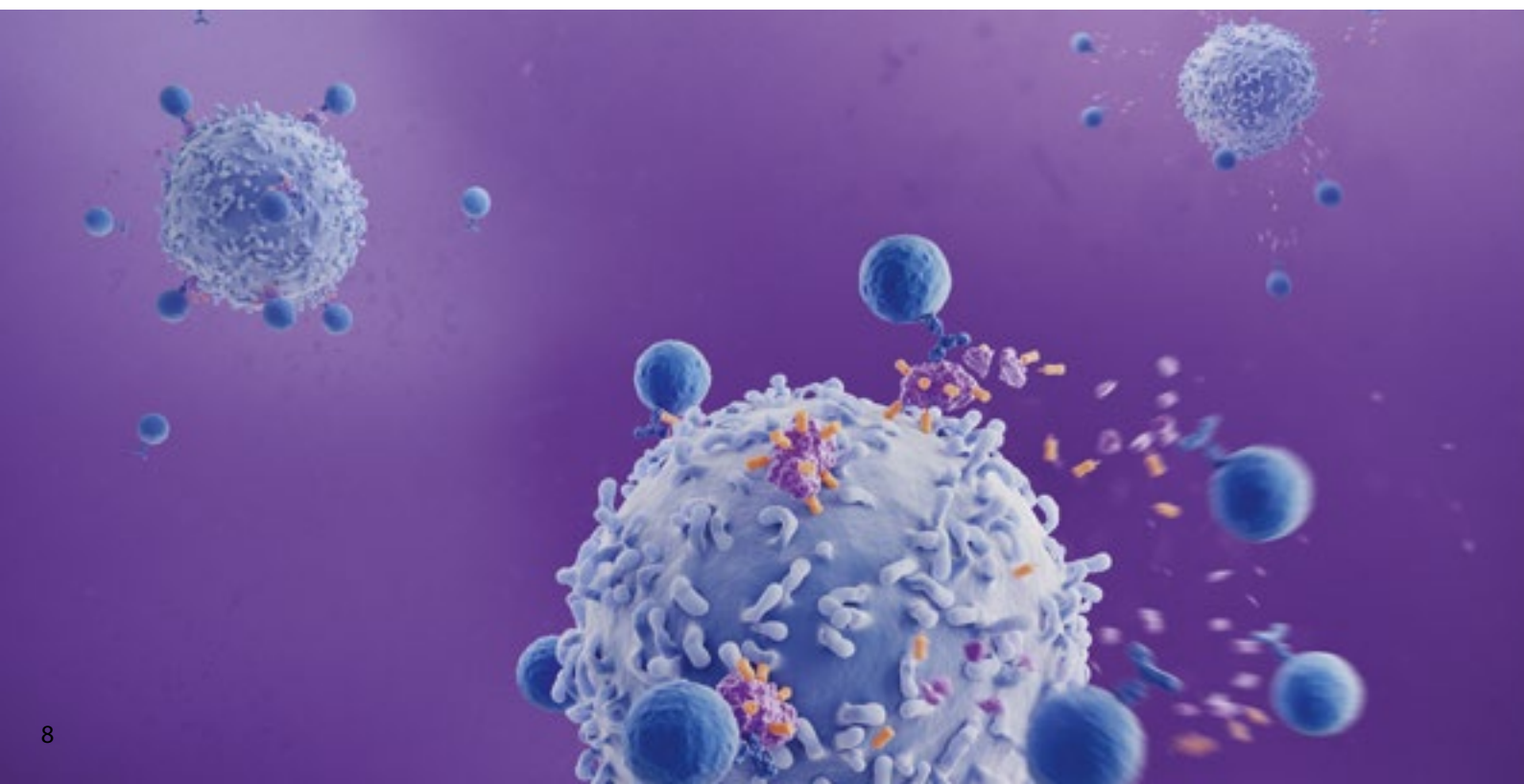
Product tables

Cell separation reagents

Products	Order no.
Anti-Mouse IgG1 MicroBeads	130-047-101
Anti-Mouse IgG2a+b MicroBeads	130-047-201
Anti-Mouse IgM MicroBeads	130-047-301
B-1a Cell Isolation Kit	130-097-413
B Cell Isolation Kit	130-090-862
CD5 (Ly-1) MicroBeads	130-049-301
CD19 MicroBeads	130-121-301
CD22 MicroBeads	130-103-776
CD23 MicroBeads	130-098-784
CD43 (Ly-48) MicroBeads	130-049-801
CD45R (B220) MicroBeads	130-049-501
CD93 MicroBeads	130-105-753
CD138 MicroBeads	130-098-257
CD138 ⁺ Plasma Cell Isolation Kit	130-092-530
Germinal Center B Cell (PNA) MicroBead Kit	130-110-479
Memory B Cell Isolation Kit	130-095-838
MZ and FO B Cell Isolation Kit	130-100-366
Pan B Cell Isolation Kit II	130-104-443
Regulatory B Cell Isolation Kit	130-095-873

Cell culture reagents

Products	Research grade	Premium grade	Functional grade
Flt3-Ligand	●		
IFN- γ	●	●	
IL-4	●	●	
IL-5	●		
IL-6	●	●	
IL-13	●		
SCF	●	●	
Stimulation antibodies			
CD2 antibody pure			●
CD48 antibody pure			●
CD127 antibody pure			●
CD134 (OX40) antibody pure			●
CD137 antibody pure			●
CD153 antibody pure			●
CD252 (OX40L) antibody pure			●
CD275 (B7-H2) antibody pure			●
CD279 (PD1) antibody pure			●



Flow cytometry antibodies

Antigen	Hybridoma Clones	REAffinity Clones	REALease Clones
CD1d	1B1		
CD5	53-7.3	REA421	REAL174
CD11b	M1/70.15.11.5	REA592	REAL113
CD19	6D5	REA749	
CD21/CD35	7E9	REA800	
CD22	Cy34.1	REA1187	
CD23	B3B4	REA1068	
CD24	M1/69	REA743	
CD25	7D4	REA568	
CD27	LG.3A10	REA499	
CD38	90.4	REA616	REAL181
CD43	L11	REA840	
CD45	30F11	REA737	REAL180
CD45.1	A20	REA1179	
CD45.2	104-2	REA1223	
CD45R (B220)	RA3-6B2	REA755	REAL132
CD45RA	14.8	REA639	
CD45RB	C363-16A	REA1180	
CD80	16-10A1	REA983	
CD86	PO3.3	REA1190	
CD93		REA298	
CD95 (FAS)		REA453	
CD138		REA104	
IL-10	JES5-16E3	REA1008	
IgD	11-26c.2a	REA772	
IgM	X-54	REA979	
IgG1	X-56	REA1017	
IgG2ab	X-57		
MHC Class II variants	M5/114.15.2	REA510	REAL362
MHC Class II variants	M5/114.15.2	REA528	REAL362
MHC Class II variants	M5/114.15.2	REA564	REAL362
MHC Class II variants	M5/114.15.2	REA610	REAL362
MHC Class II variants	M5/114.15.2	REA813	REAL362

Flow cytometry cocktails and kits

Products	Order no.
Mouse IL-2 Secretion Assay – Detection Kit (APC)	130-090-987
Mouse IL-2 Secretion Assay – Detection Kit (PE)	130-090-491
Mouse IL-4 Secretion Assay – Detection Kit (PE)	130-090-479
Mouse IL-5 Secretion Assay – Detection Kit (APC)	130-091-174
Mouse IL-5 Secretion Assay – Detection Kit (PE)	130-091-166
Mouse IL-10 Secretion Assay – Detection Kit (APC)	130-090-939
Mouse IL-10 Secretion Assay – Detection Kit (PE)	130-090-489
Mouse IFN- γ Secretion Assay – Detection Kit (APC)	130-090-984
Mouse IFN- γ Secretion Assay – Detection Kit (PE)	130-090-516

MACSPlex Custom Cytokine Assays

Products	Order no.
MACSPlex Cytokine Basic Kit, human and mouse	130-109-701
MACSPlex Cytokine 10 Kit	130-101-740
MACSPlex GM-CSF Reagents Kit	130-109-702
MACSPlex IFN- γ Reagents Kit	130-109-687
MACSPlex IL-2 Reagents Kit	130-109-686
MACSPlex IL-4 Reagents Kit	130-109-690
MACSPlex IL-5 Reagents Kit	130-109-692
MACSPlex IL-10 Reagents Kit	130-109-698
MACSPlex IL-12 Reagents Kit	130-109-696
MACSPlex IL-17A Reagents Kit	130-109-685
MACSPlex IL-23 Reagents Kit	130-109-700
MACSPlex TNF- α Reagents Kit	130-109-703

For more information about antibodies and fluorescent conjugates for flow cytometry visit:
www.miltenyibiotec.com/flowcytometry

Other species

Product tables

Cell separation reagents

Products	Order no.
Rat	
Anti-Rat Kappa MicroBeads	130-047-401
CD43 MicroBeads	130-090-742
CD45R MicroBeads	130-090-495
Non-human primate	
CD20 MicroBeads	130-091-105

Flow cytometry antibodies

Antigen	Hybridoma Clones	REAffinity Clones
Rat		
CD5		REA750
CD24		REA405
CD32		REA256
CD38		REA683
CD43		REA503
CD45		REA504
CD45R		REA450
CD45RC		REA1000
CD86	24F	
CD200		REA532
CD252	ATM-2	
IgG1	RG11/39.2	
IgG2a	RG7/1.30	
IgG2b	RG7/11.1	
MHC Class II		REA510
Non-human primate		
CD45	MB4-6D6	REA1023

Visit our webpage for our complete product portfolios and scientific resources

- Webinars
- Videos
- Protocols and workflows
- Application notes
- Posters
- and more ...

► **[miltenyibiotec.com](https://www.miltenyibiotec.com)**





Miltenyi Biotec

Germany/Austria

Miltenyi Biotec B.V. & Co. KG
Friedrich-Ebert-Straße 68
51429 Bergisch Gladbach
Germany
Phone +49 2204 8306-0
Fax +49 2204 85197
macsde@miltenyi.com

USA/Canada

Miltenyi Biotec, Inc.
2303 Lindbergh Street
Auburn, CA 95602, USA
Phone 800 FOR MACS
Phone +1 866 811 4466
Fax +1 877 591 1060
macsau@miltenyi.com

Australia

Miltenyi Biotec
Australia Pty. Ltd.
Unit 11, 2 Eden Park Drive
Macquarie Park, NSW 2113
Australia
Phone +61 2 8877 7400
Fax +61 2 9889 5044
macsau@miltenyi.com

Benelux

Miltenyi Biotec B.V.
Sandifordreef 17
2333 ZZ Leiden
The Netherlands
macsnl@miltenyi.com
Customer service

The Netherlands
Phone 0800 4020120
Fax 0800 4020100

Customer service Belgium

Phone 0800 94016
Fax 0800 99626

Customer service Luxembourg

Phone 800 24971
Fax 800 24984

China

Miltenyi Biotec Technology &
Trading (Shanghai) Co., Ltd.
Room A401-402, 4/F
No. 1077, Zhangheng Road
Pudong New Area
201203 Shanghai, P.R. China
Phone +86 21 6235 1005-0
Fax +86 21 6235 0953
macscn@miltenyi.com

France

Miltenyi Biotec SAS
10 rue Mercœur
75011 Paris, France
Phone +33 1 56 98 16 16
macsfr@miltenyi.com

Hong Kong

Miltenyi Biotec Hong Kong Ltd.
Unit 301, Lakeside 1
No. 8 Science Park West Avenue
Hong Kong Science Park
Pak Shek Kok, New Territories
Hong Kong
Phone +852 3751 6698
Fax +852 3619 5772
macshk@miltenyi.com

Italy

Miltenyi Biotec S.r.l.
Via Paolo Nanni Costa, 30
40133 Bologna
Italy
Phone +39 051 6 460 411
Fax +39 051 6 460 499
macsit@miltenyi.com

Japan

Miltenyi Biotec K.K.
NEX-Eitai Building 5F
16-10 Fuyuki, Koto-ku
Tokyo 135-0041, Japan
Phone +81 3 5646 8910
Fax +81 3 5646 8911
macsjp@miltenyi.com

Nordics and Baltics

Miltenyi Biotec Norden AB
Medicon Village
Scheeleorget 1
223 81 Lund
Sweden
macsse@miltenyi.com
Customer service Sweden
Phone 0200 111 800
Fax +46 280 72 99

Customer service Denmark

Phone 80 20 30 10
Fax +46 46 280 72 99

Customer service

**Norway, Finland, Iceland,
and Baltic countries**
Phone +46 46 280 72 80
Fax +46 46 280 72 99

Singapore

Miltenyi Biotec Asia Pacific Pte Ltd.
438B Alexandra Road, Block B
Alexandra Technopark
#06-01
Singapore 119968
Phone +65 6238 8183
Fax +65 6238 0302
macssg@miltenyi.com

South Korea

Miltenyi Biotec Korea Co., Ltd.
Arigi Bldg. 8F
562 Nonhyeon-ro
Gangnam-gu
Seoul 06136, South Korea
Phone +82 2 555 1988
Fax +82 2 555 8890
macskr@miltenyi.com

Spain

Miltenyi Biotec S.L.
C/Luis Buñuel 2
Ciudad de la Imagen
28223 Pozuelo de Alarcón (Madrid)
Spain
Phone +34 91 512 12 90
Fax +34 91 512 12 91
macses@miltenyi.com

Switzerland

Miltenyi Biotec Swiss AG
Gibelinstrasse 27
4500 Solothurn
Switzerland
Phone +41 32 623 08 47
Fax +49 2204 85197
macsch@miltenyi.com

United Kingdom

Miltenyi Biotec Ltd.
Almac House, Church Lane
Bisley, Surrey GU24 9DR, UK
Phone +44 1483 799 800
Fax +44 1483 799 811
macsuk@miltenyi.com

www.miltenyibiotec.com

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

This publication is provided for information purposes only, and the contents hereof are subject to change without notice. While Miltenyi Biotec has taken great care and attention to ensure the quality of the information contained herein at the time of publication, no warranties are provided that the content of this publication is accurate, complete, reliable, up to date, and error-free. Miltenyi Biotec assumes no responsibility or liability arising out of the application or use of any information described herein except as expressly confirmed in writing by authorized representative of Miltenyi Biotec. Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. MACS, MACSxpress, the Miltenyi Biotec logo, PepTivator, REAfinity, REAlease, and StraightFrom are registered trademarks or trademarks of Miltenyi Biotec and/or its affiliates in various countries worldwide. Leukopak is a registered trademark of StemExpress LLC. Copyright © 2022 Miltenyi Biotec and/or its affiliates.