

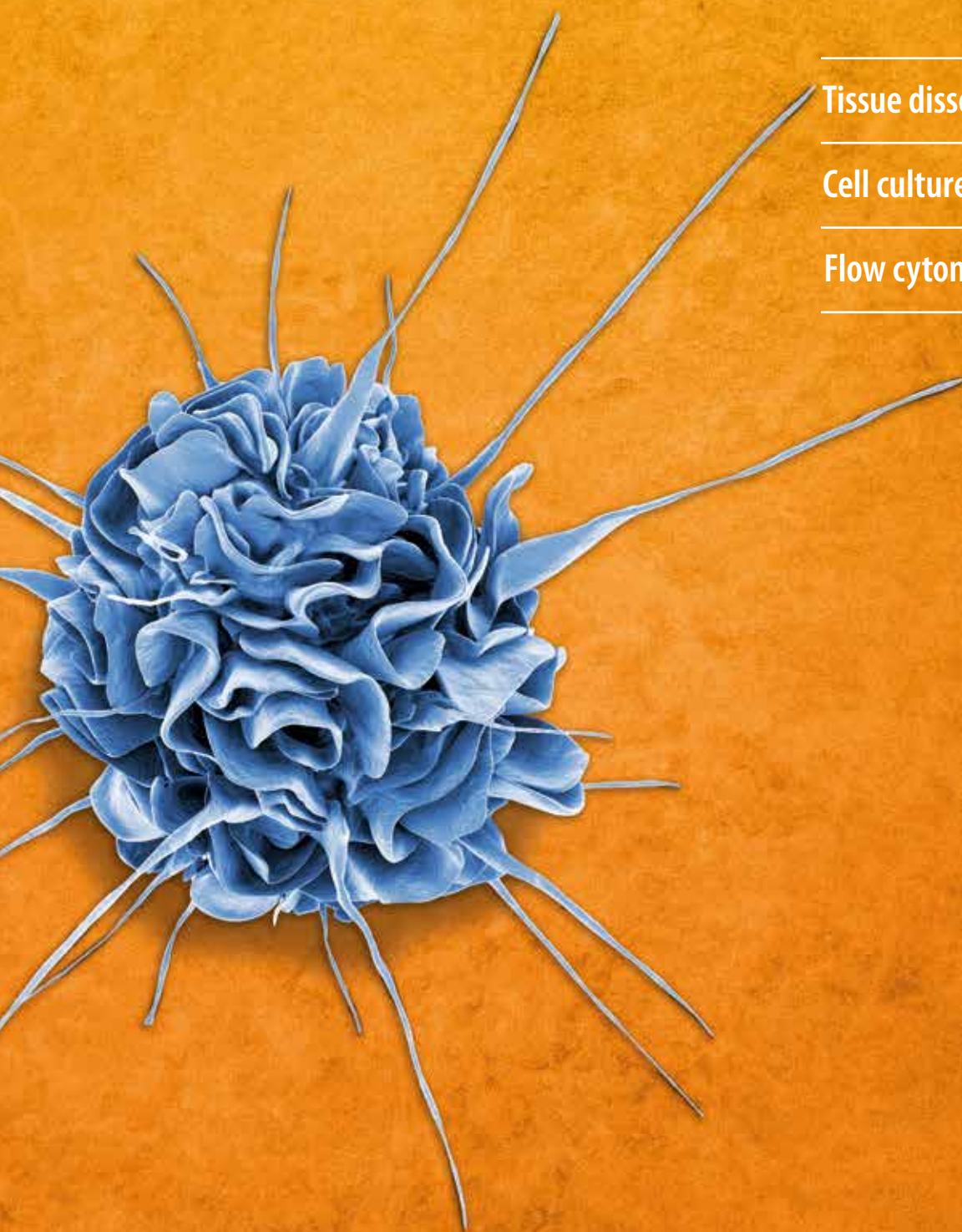
Product quick guide

All you need for myeloid cell research

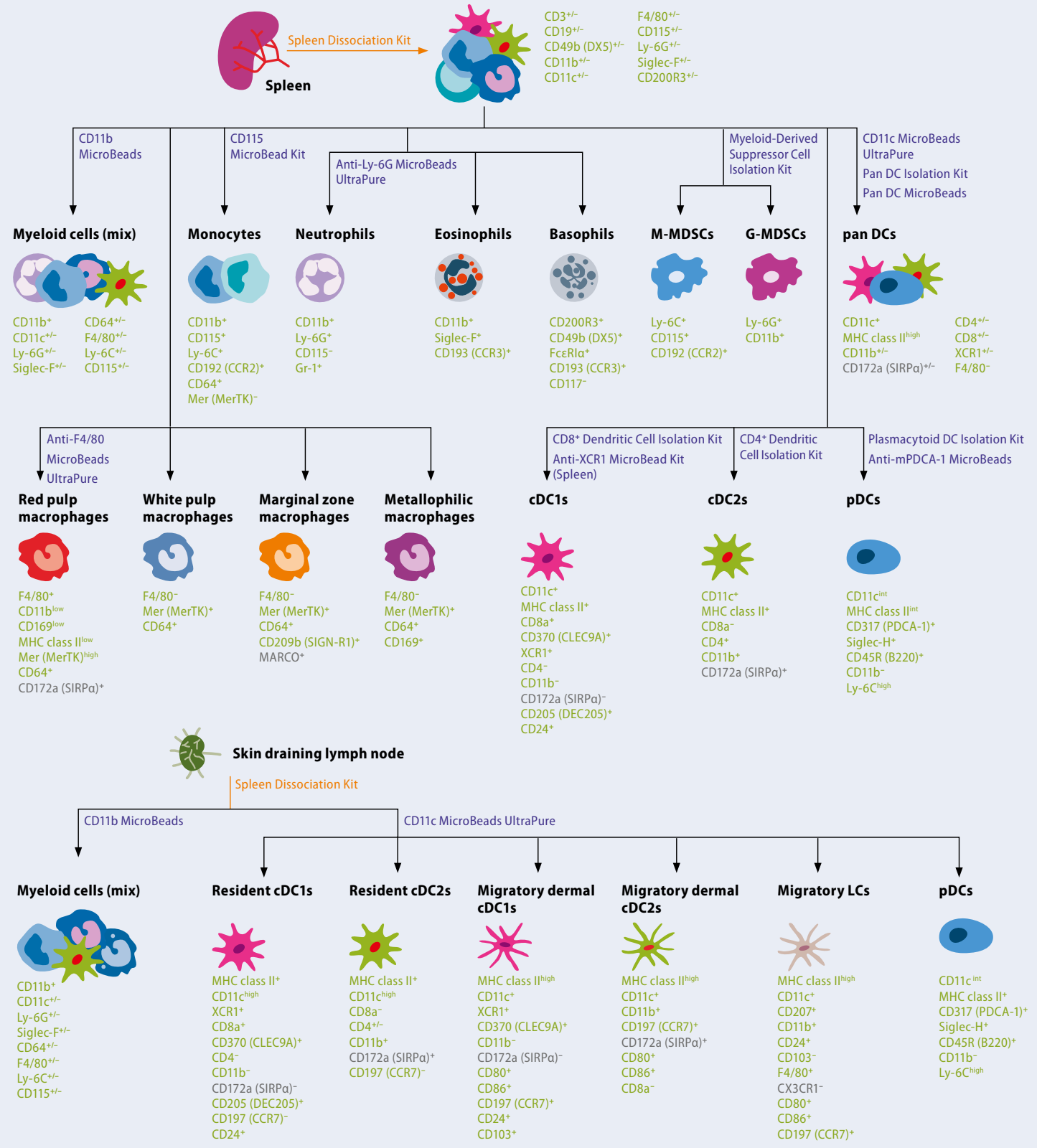
Tissue dissociation & cell separation

Cell culture & stimulation

Flow cytometry



Mouse myeloid cells from spleen and lymph nodes



Legend

→ separation
→ dissociation

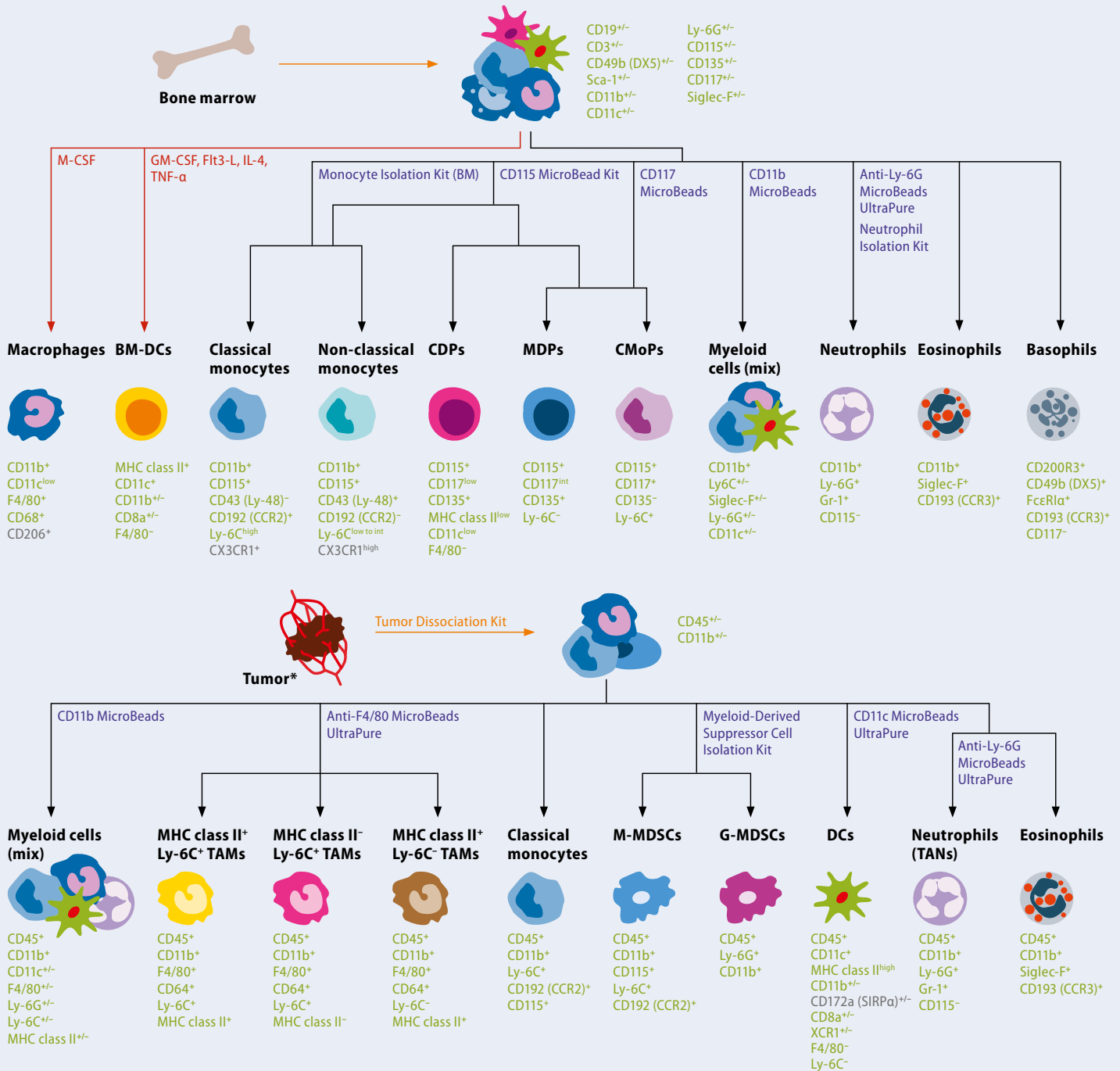
- MACS[®] Sample Preparation products (RUO)
- MACS Cell Separation products (RUO)
- MACS Cell Culture and Stimulation products (RUO)
- MACS Flow Cytometry products (RUO)

DC: dendritic cell
cDC: conventional or myeloid DC
pDC: plasmacytoid DC
M-MDSC: monocytic myeloid-derived suppressor cell
G-MDSC: granulocytic myeloid-derived suppressor cell
LC: Langerhans cells

- no expression
^{low} low expression
^{int} intermediate expression
⁺ expression

^{high} high expression
⁺ expression dependent on subset

Mouse myeloid cells from bone marrow and tumor



Legend

- separation
- differentiation
- dissociation

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

- DC: dendritic cell
- cDC: conventional or myeloid DC
- pDC: plasmacytoid DC
- BM: bone marrow
- BM-DC: bone marrow-derived DC
- CDP: common dendritic cell progenitor
- MDP: monocyte and DC progenitor

- CMoP: common monocyte progenitor
- M-MDSC: monocytic myeloid-derived suppressor cell
- G-MDSC: granulocytic myeloid-derived suppressor cell
- LC: Langerhans cell
- TAM: tumor-associated macrophage
- TAN: tumor-associated neutrophil

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

* Please contact our technical support to know which tumor types were tested with indicated products.

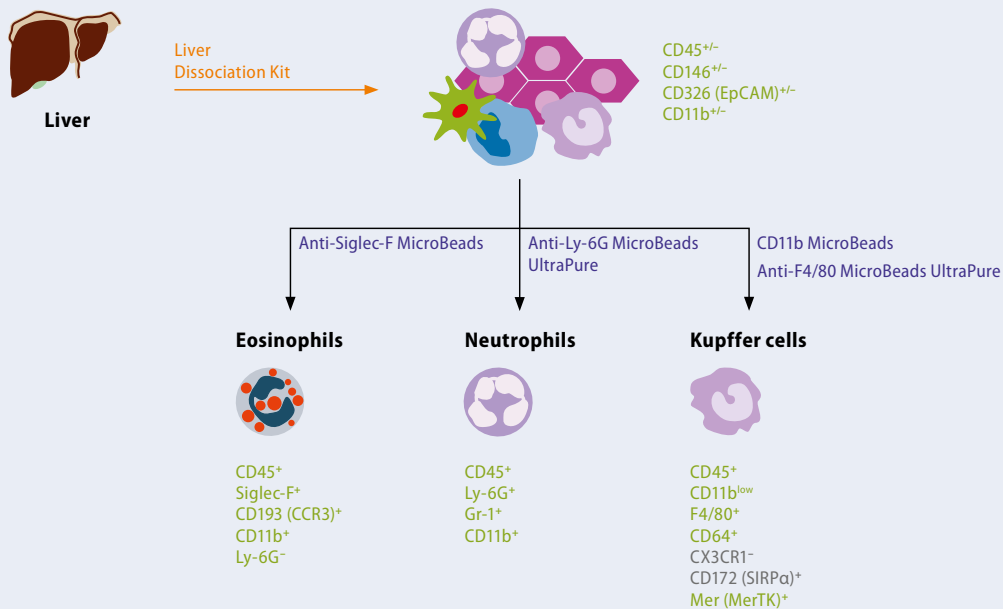
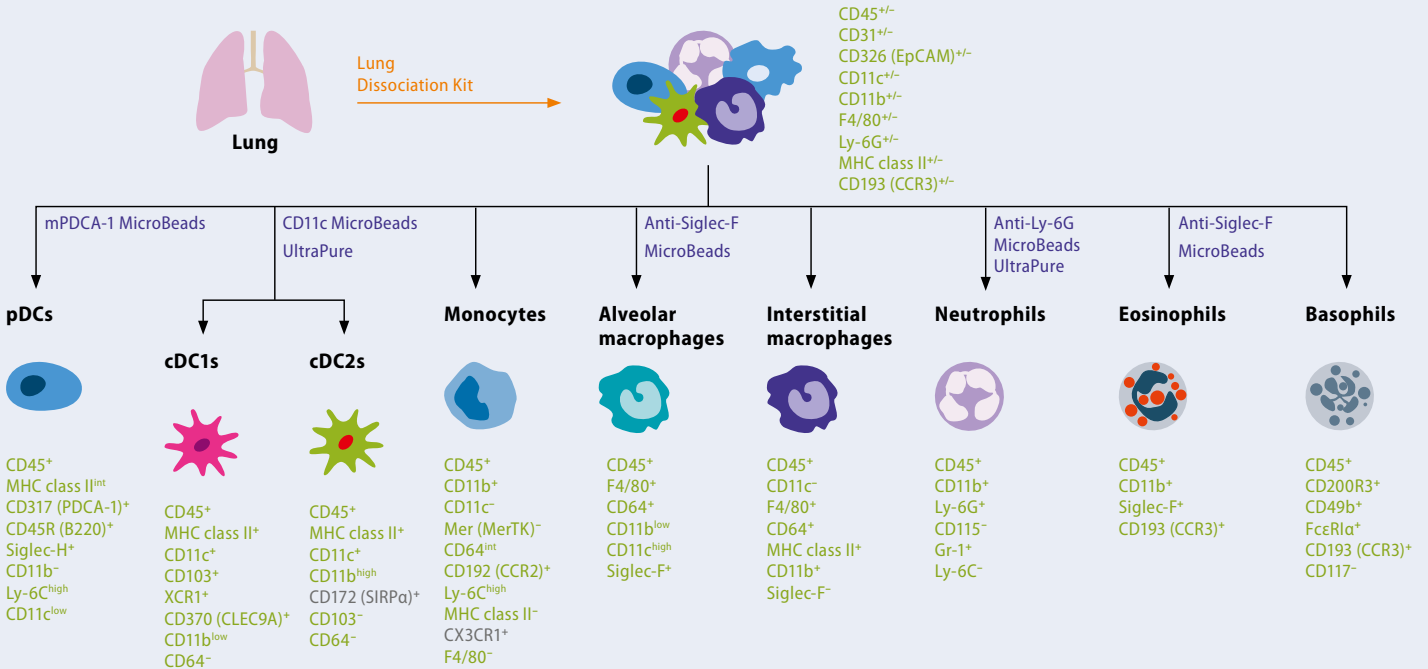
Miltenyi Biotec B.V. & Co. KG | Friedrich-Ebert-Straße 68 | 51429 Bergisch Gladbach | Germany | Phone +49 2204 8306-0 | Fax +49 2204 85197
 macs@miltenyibiotec.de | www.miltenyibiotec.com

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

[miltenyibiotec.com](http://www.miltenyibiotec.com)

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, MACSexpress, StraightFrom, and the MACS logo are registered trademarks or trademarks of Miltenyi Biotec B.V. & Co. KG and/or its affiliates in various countries worldwide. All other trademarks mentioned in this document are the property of their respective owners and are used for identification purposes only. Leukopak is a registered trademark of StemExpress, LCC. Copyright © 2019 Miltenyi Biotec B.V. & Co. KG and/or its affiliates. All rights reserved.

Mouse myeloid cells from lung and liver



Legend

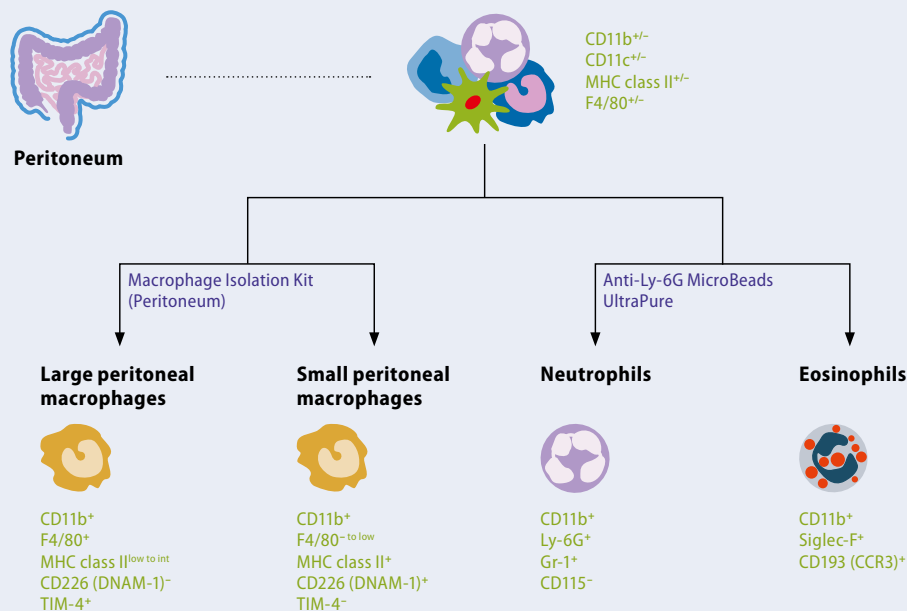
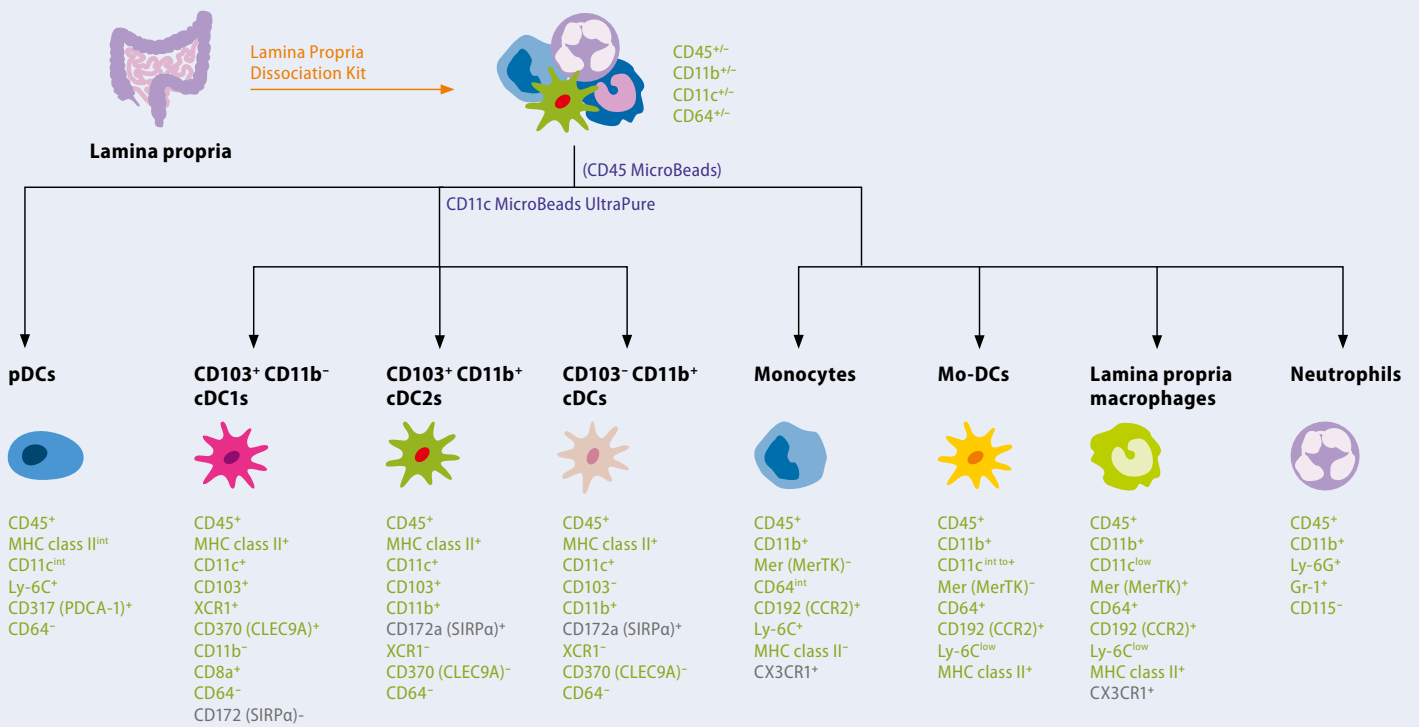
→ separation
→ dissociation

DC: dendritic cell
cDC: conventional or myeloid DC
pDC: plasmacytoid DC

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

■ MACS[®] Sample Preparation products (RUO)
■ MACS Cell Separation products (RUO)
■ MACS Cell Culture and Stimulation products (RUO)
■ MACS Flow Cytometry products (RUO)

Mouse myeloid cells from lamina propria and peritoneum



Legend

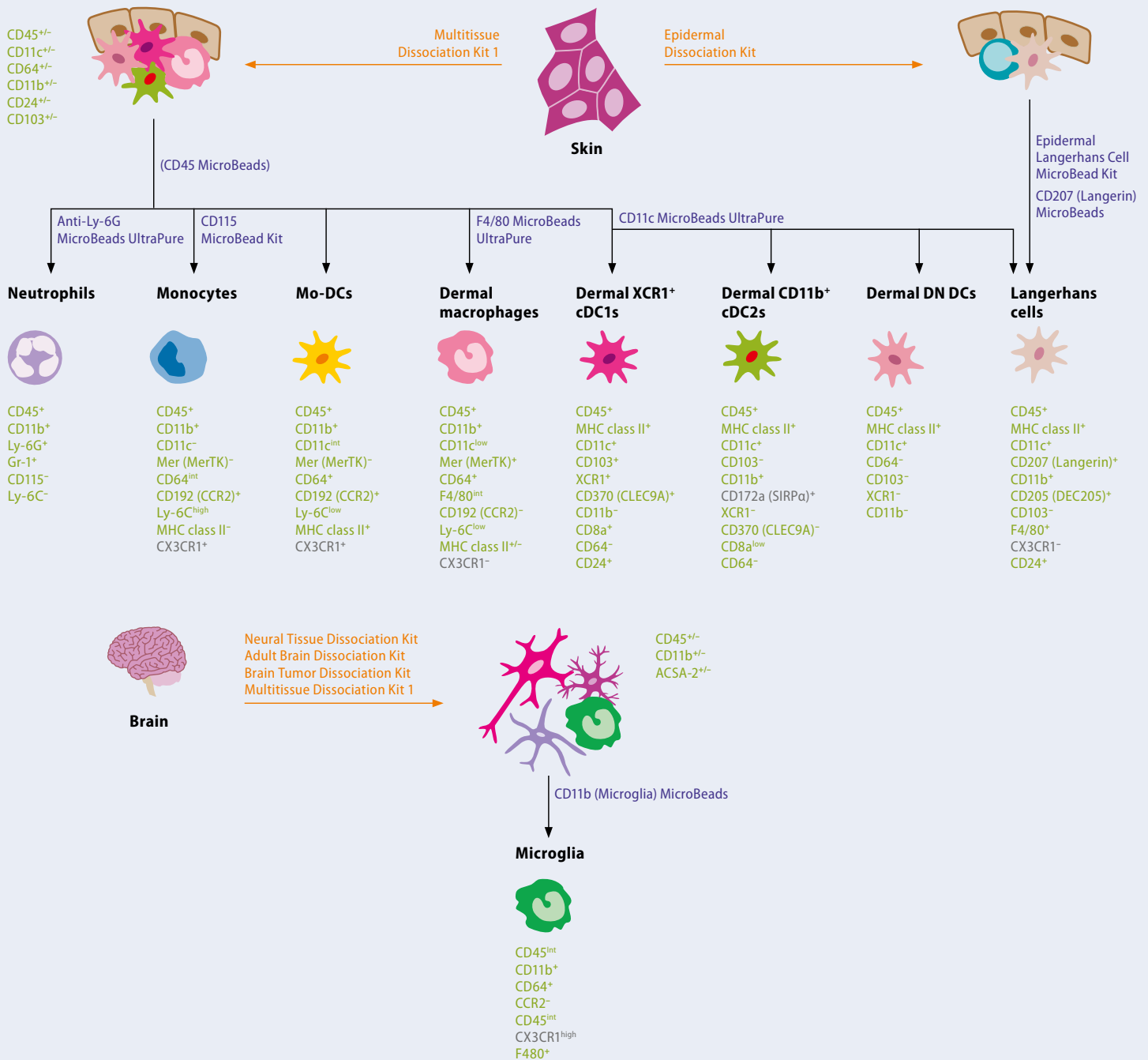
→ separation
 → dissociation

■ MACS[®] Sample Preparation products (RUO)
 ■ MACS Cell Separation products (RUO)
 ■ MACS Cell Culture and Stimulation products (RUO)
 ■ MACS Flow Cytometry products (RUO)

DC: dendritic cell
cDC: conventional or myeloid DC
pDC: plasmacytoid DC
Mo-DC: monocyte-derived DC

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

Mouse myeloid cells from skin and brain



Legend

- separation
- dissociation

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

DC: dendritic cell
cDC: conventional or myeloid DC
Mo-DC: monocyte derived DC
Dermal DN: Dermal double negative

- no expression
^{low} low expression
^{int} intermediate expression
⁺ expression

^{high} high expression
^{+/-} expression dependent on subset

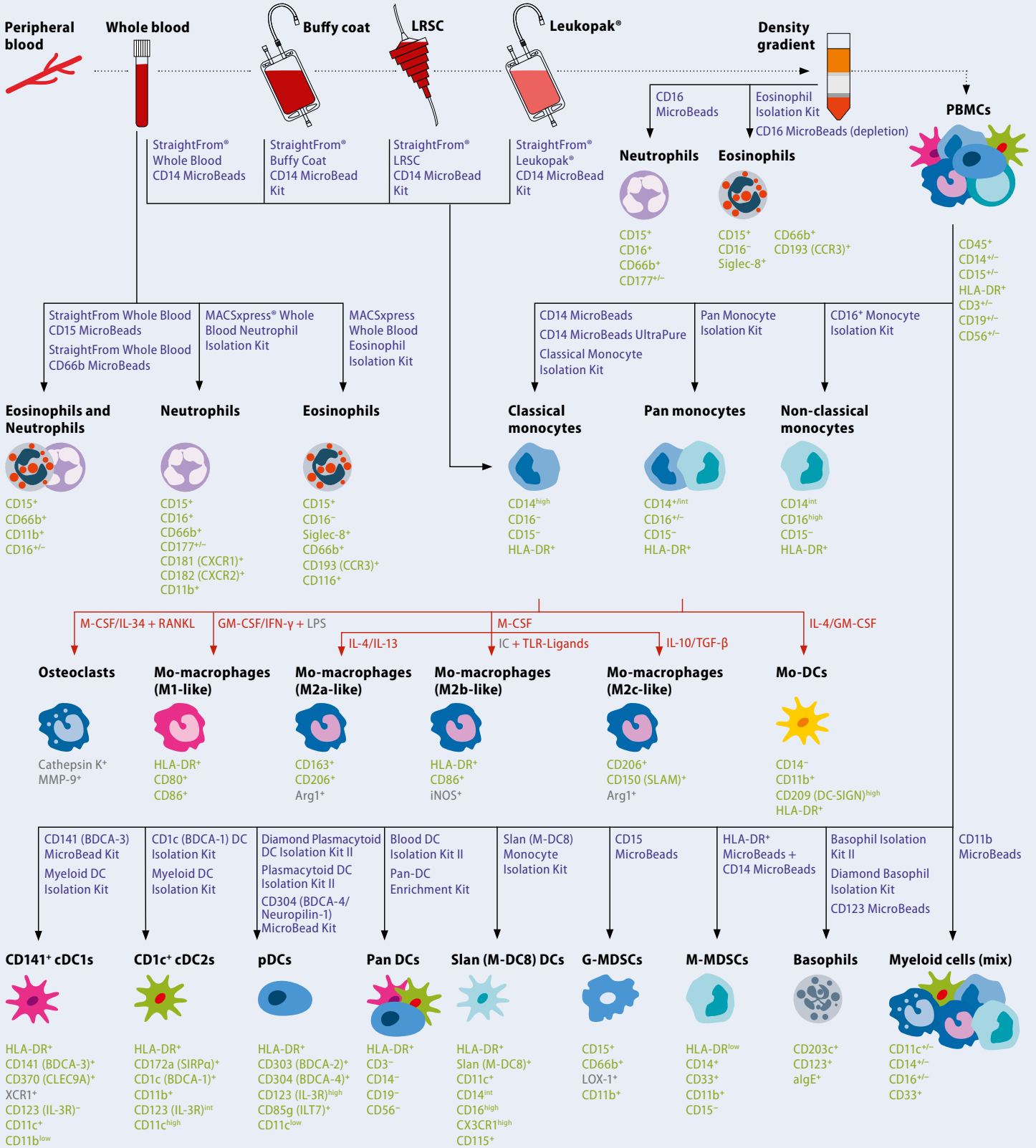
Miltenyi Biotec B.V. & Co. KG | Friedrich-Ebert-Straße 68 | 51429 Bergisch Gladbach | Germany | Phone +49 2204 8306-0 | Fax +49 2204 85197
 macs@miltenyibiotec.de | www.miltenyibiotec.com

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

▶ miltenyibiotec.com

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, StraightFrom, and the MACS logo are registered trademarks or trademarks of Miltenyi Biotec B.V. & Co. KG and/or its affiliates in various countries worldwide. All other trademarks mentioned in this document are the property of their respective owners and are used for identification purposes only. Leukopak is a registered trademark of StemExpress, LLC. Copyright © 2019 Miltenyi Biotec B.V. & Co. KG and/or its affiliates. All rights reserved.

Human myeloid cells from peripheral blood



Legend

- separation
- differentiation (cultivation)

- MACS® Cell Separation products (RUO)
- MACS Cell Culture and Stimulation products (RUO)
- MACS Flow Cytometry products (RUO)

- DC:** dendritic cell
- cDC:** conventional or myeloid DC
- pDC:** plasmacytoid DC
- Mo-DC:** monocyte-derived DC
- PBMC:** peripheral blood mononuclear cell

- M-MDSC:** monocytic myeloid-derived suppressor cell
- G-MDSC:** granulocytic myeloid-derived suppressor cell
- IC:** immunocomplexes

- no expression
- low low expression
- int intermediate expression
- + expression

- ^{high} high expression
- ^{+/-} expression dependent on subset



Miltenyi Biotec

Germany/Austria

Miltenyi Biotec B.V. & Co. KG
Friedrich-Ebert-Str. 68
51429 Bergisch Gladbach
Germany
Phone +49 2204 8306-0
Fax +49 2204 85197
macs@miltenyibiotec.de

USA/Canada

Miltenyi Biotec Inc.
2303 Lindbergh Street
Auburn, CA 95602, USA
Phone 800 FOR MACS
Phone +1 530 888 8871
Fax +1 877 591 1060
macs@miltenyibiotec.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
macs@miltenyibiotec.com.au

Benelux

Miltenyi Biotec B.V.
Sandifortdreef 17
2333 ZZ Leiden
The Netherlands
macs@miltenyibiotec.nl

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 401
No. 1077, Zhangheng Road
Pudong New Area
201203 Shanghai, P.R. China
Phone +86 21 6235 1005
Fax +86 21 6235 0953
macs@miltenyibiotec.com.cn

France

Miltenyi Biotec SAS
10 rue Mercoeur
75011 Paris, France
Phone +33 1 56 98 16 16
Fax +33 1 56 98 16 17
macs@miltenyibiotec.fr

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
macs@miltenyibiotec.it

Japan

Miltenyi Biotec K.K.
Nittsu-Eitai Building 5F
16-10 Fuyuki, Koto-ku
Tokyo 135-0041, Japan
Phone +81 3 5646 8910
Fax +81 3 5646 8911
macs@miltenyibiotec.jp

Nordics and Baltics

Miltenyi Biotec Norden AB
Scheelevägen 17
223 70 Lund
Sweden
macs@miltenyibiotec.se

Customer service Sweden

Phone 0200 111 800
Fax 046 280 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.
100 Beach Road
#28-06 to 28-08 Shaw Tower
Singapore 189702
Phone +65 6238 8183
Fax +65 6238 0302
macs@miltenyibiotec.com.sg

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
macs@miltenyibiotec.co.kr

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
macs@miltenyibiotec.es

Switzerland

Miltenyi Biotec Swiss AG
Gibelinstrasse 27
4500 Solothurn
Switzerland
Phone +41 32 623 08 47
Fax +49 2204 85197
macs@miltenyibiotec.ch

United Kingdom

Miltenyi Biotec Ltd.
Almac House, Church Lane
Bisley, Surrey GU24 9DR, UK
Phone +44 1483 799 800
Fax +44 1483 799 811
macs@miltenyibiotec.co.uk

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, StraightFrom, and the MACS logo are registered trademarks or trademarks of Miltenyi Biotec B.V. & Co. KG and/or its affiliates. All other trademarks mentioned in this document are the property of their respective owners and are used for identification purposes only. Leukopak is a registered trademark of StemExpress, LCC. Copyright © 2019 Miltenyi Biotec B.V. & Co. KG and/or its affiliates. All rights reserved.