

# The broadest portfolio for your dendritic cell research

---

Tissue dissociation and cell separation

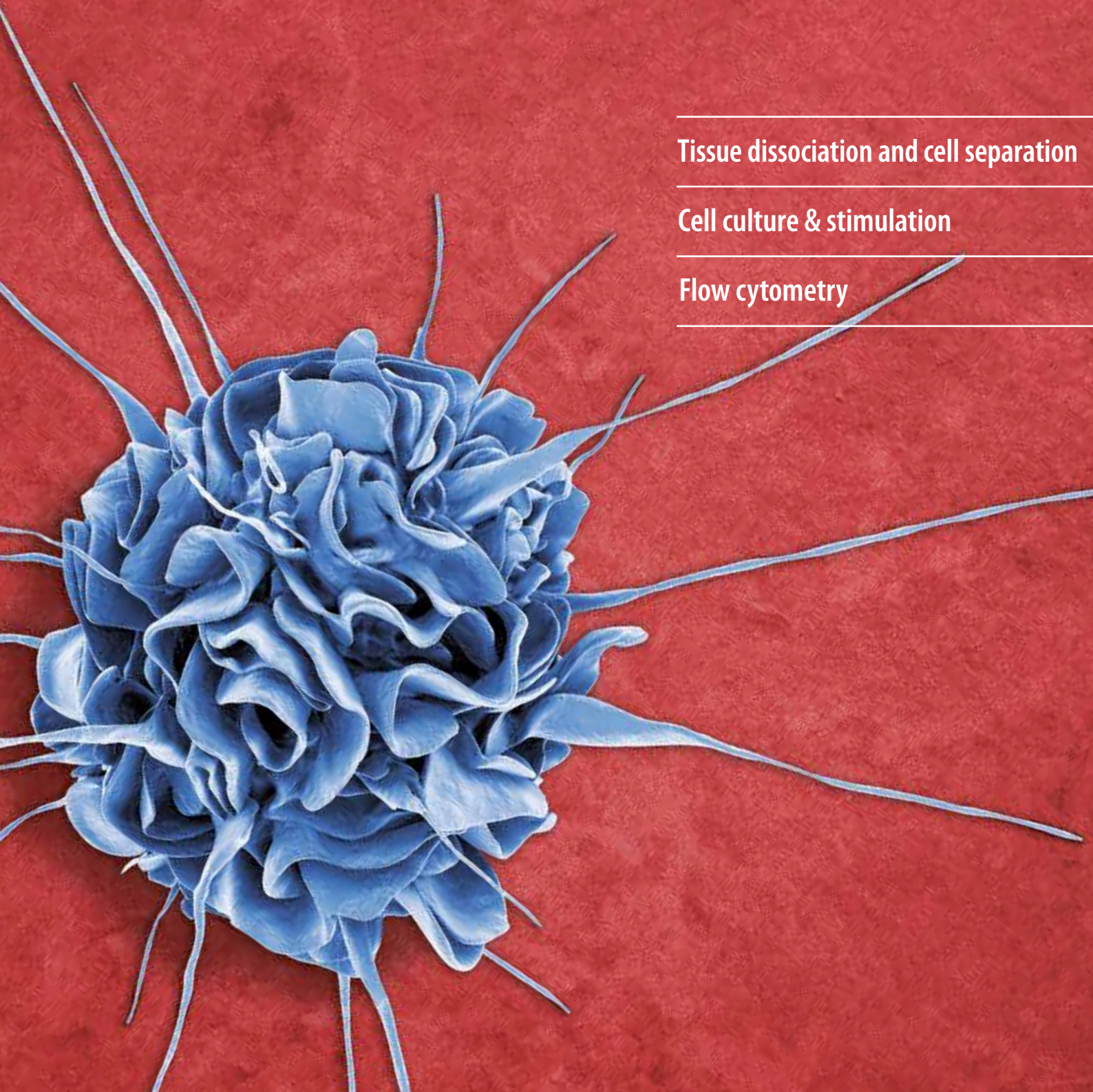
---

Cell culture & stimulation

---

Flow cytometry

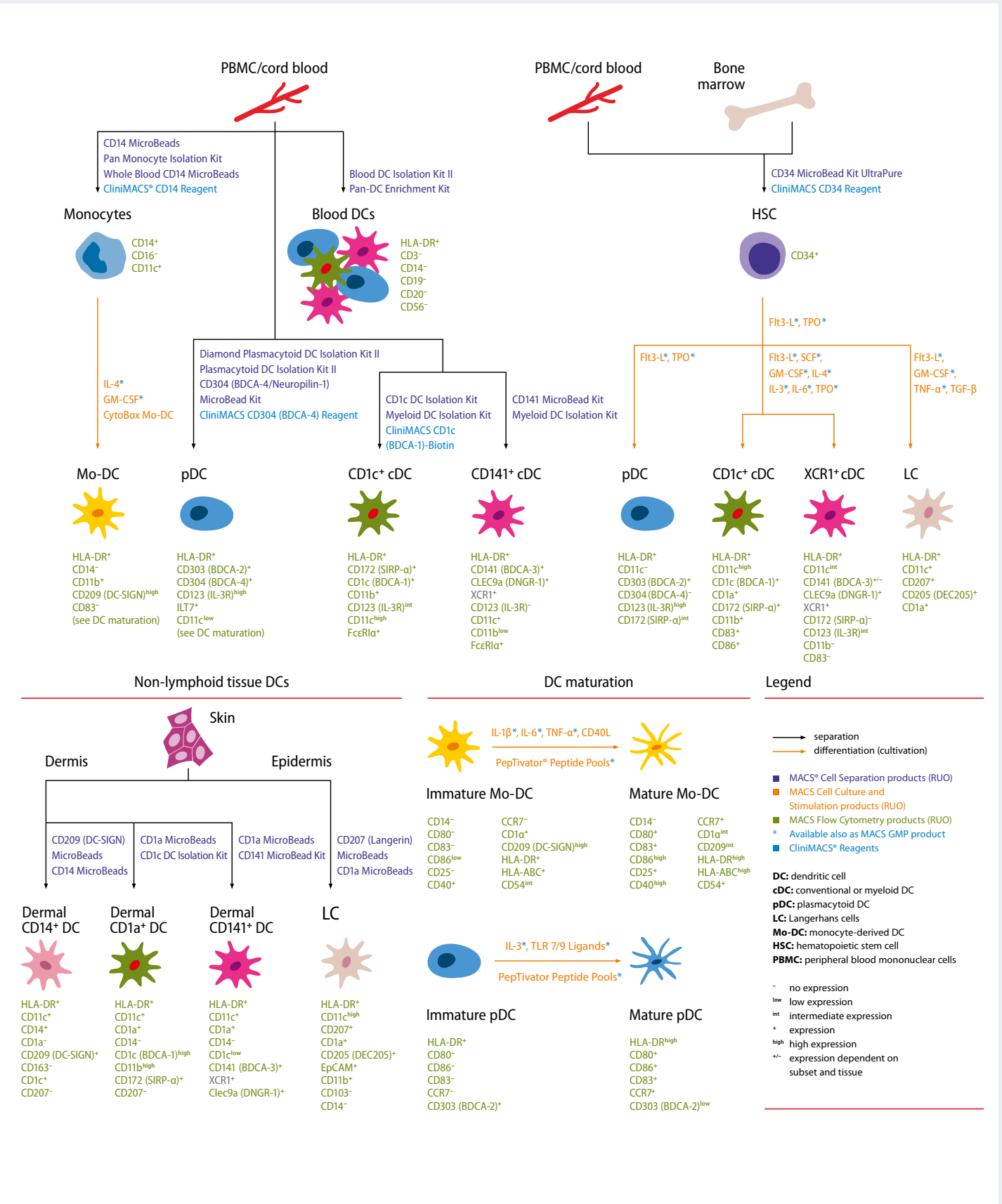
---





# Human dendritic cell generation

## Overview



# Human

## Product tables

### Cell separation reagents

Products	Order no.
CD14 MicroBeads	130-050-201
Pan Monocyte Isolation Kit	130-096-537
Whole Blood CD14 MicroBeads	130-090-879
Blood Dendritic Cell Isolation Kit II	130-091-379
Pan-DC Enrichment Kit	130-100-777
Diamond Plasmacytoid Dendritic Cell Isolation Kit II	130-097-240
Plasmacytoid Dendritic Cell Isolation Kit II	130-097-415
CD304 (BDCA-4/Neuropilin-1) MicroBead Kit	130-090-532
CD1c (BDCA-1)+ Dendritic Cell Isolation Kit	130-090-506
Myeloid Dendritic Cells Isolation Kit	130-094-487
CD141 (BDCA-3) MicroBead Kit	130-090-512
CD207 (Langerin) MicroBeads	130-097-898
CD1a MicroBeads	130-051-001
CD209 (DC-SIGN) MicroBead Kit	130-092-868

For more information visit:

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

### Cell culture reagents

Products	Research grade	Premium grade	MACS® GMP
Human IL-4	●	●	●
Human GM-CSF	●	●	●
Human IL-6	●	●	●
Human IL-1β	●	●	●
Human TNF-α	●	●	●
CytoBox Mo-DC		●	
Human SCF	●	●	●
Human Flt3-Ligand	●	●	●
Human TPO	●	●	●
Human IL-3	●	●	●
Human TGF-β1		●	
TLR9-Ligands	●		●
TLR7/8-Ligands	●		
PepTivators (Peptide pools)	●	●	●

For more information about custom fillings and cocktails visit:

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

### Flow cytometry antibodies

Antigen	Vio-Blue®	Vio-Green™	FITC	VioBright™ FITC	PE	PerCP	PerCP-Vio® 700	PE-Vio 615	PE-Vio 770	APC	APC-Vio 770	Biotin	REAlity
<b>Mo/Mo-DC</b>													
CD14		●	●		●	●	●	●	●	●	●	●	
CD16	●	●	●		●		●	●	●	●	●	●	●
CD11c	●		●	●	●		●		●	●	●	●	●
HLA-DR	●	●	●		●	●	●		●	●	●	●	●
HLA-ABC	●	●	●		●		●	●	●	●	●	●	●
CD80					●				●	●		●	●
CD86	●		●		●		●		●	●		●	●
CD83					●				●	●		●	●
CCR7	●		●		●		●		●	●		●	●
CD25				●	●					●		●	●
CD40					●				●	●		●	●
CD54					●			●	●	●		●	●
CD1a	●		●		●				●	●	●	●	●
CD209 (DC-SIGN)	●		●		●		●		●	●	●	●	●
<b>Pan DC</b>													
HLA-DR	●	●	●		●	●	●	●	●	●	●	●	●
CD3	●	●	●		●	●	●	●	●	●	●	●	●
CD14	●	●	●		●	●	●	●	●	●	●	●	●
CD19	●	●	●	●	●		●	●	●	●	●	●	●
CD56			●	●	●		●	●	●	●	●	●	●
CD20	●	●	●		●			●	●	●	●	●	●

# Human

## Product tables

### Flow cytometry antibodies

Antigen	Vio-Blue®	Vio-Green™	FITC	VioBright™ FITC	PE	PerCP	PerCP-Vio® 700	PE-Vio 615	PE-Vio 770	APC	APC-Vio 770	Biotin	REAfinity
<b>pDC</b>													
HLA-DR	●	●	●		●	●	●	●	●	●	●	●	●
CD11c	●		●	●	●		●		●	●	●	●	●
CD303 (BDCA-2)	●		●	●	●		●	●	●	●	●	●	●
CD304 (BDCA-4)				●	●				●	●		●	
CD123 (IL-3R)	●	●	●		●		●	●	●	●	●	●	
CD85g (ILT7)			●		●				●	●		●	●
CD172a (SIRP-α)			●		●				●	●	●	●	●
<b>myeloid DC</b>													
HLA-DR	●	●	●		●	●	●	●	●	●	●	●	●
CD11c	●		●	●	●		●		●	●	●	●	●
CD1c (BDCA-1)			●	●	●				●	●		●	●
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
CD123 (IL-3r)	●	●	●		●		●	●	●	●	●	●	
CD141 (BDCA-3)	●		●	●	●		●		●	●	●	●	●
CLEC9A	●		●	●	●					●		●	
CD1a	●		●		●				●	●	●	●	
CD14	●	●	●		●	●	●	●	●	●	●	●	●
CD207	●		●		●				●	●	●	●	
CD172a (SIRP-α)			●		●				●	●	●	●	●
CD209 (DC-SIGN)	●		●		●		●		●	●	●	●	●
CD163	●		●		●				●	●		●	
FcεR1α			●		●		●		●	●	●		●
<b>LC</b>													
HLA-DR	●	●	●		●	●	●	●	●	●	●	●	●
CD11c	●		●	●	●		●		●	●	●	●	●
CD207	●		●		●				●	●	●	●	
CD1a	●		●		●		●		●	●	●	●	●
CD205 (DEC205)				●	●				●	●		●	
CD326 (EpCAM)	●		●		●			●	●	●	●	●	
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
CD103			●		●		●		●	●	●	●	●
CD14	●	●	●		●	●	●	●	●	●	●	●	●

For more information about MACS Antibodies and recombinantly generated REAfinity Antibodies visit [www.miltenyibiotec.com/antibodies](http://www.miltenyibiotec.com/antibodies)  
For customized fillings or bulk requests of pure antibodies for mass cytometry please go to: [www.miltenyibiotec.com/CytoF](http://www.miltenyibiotec.com/CytoF)

### Flow cytometry cocktails and kits

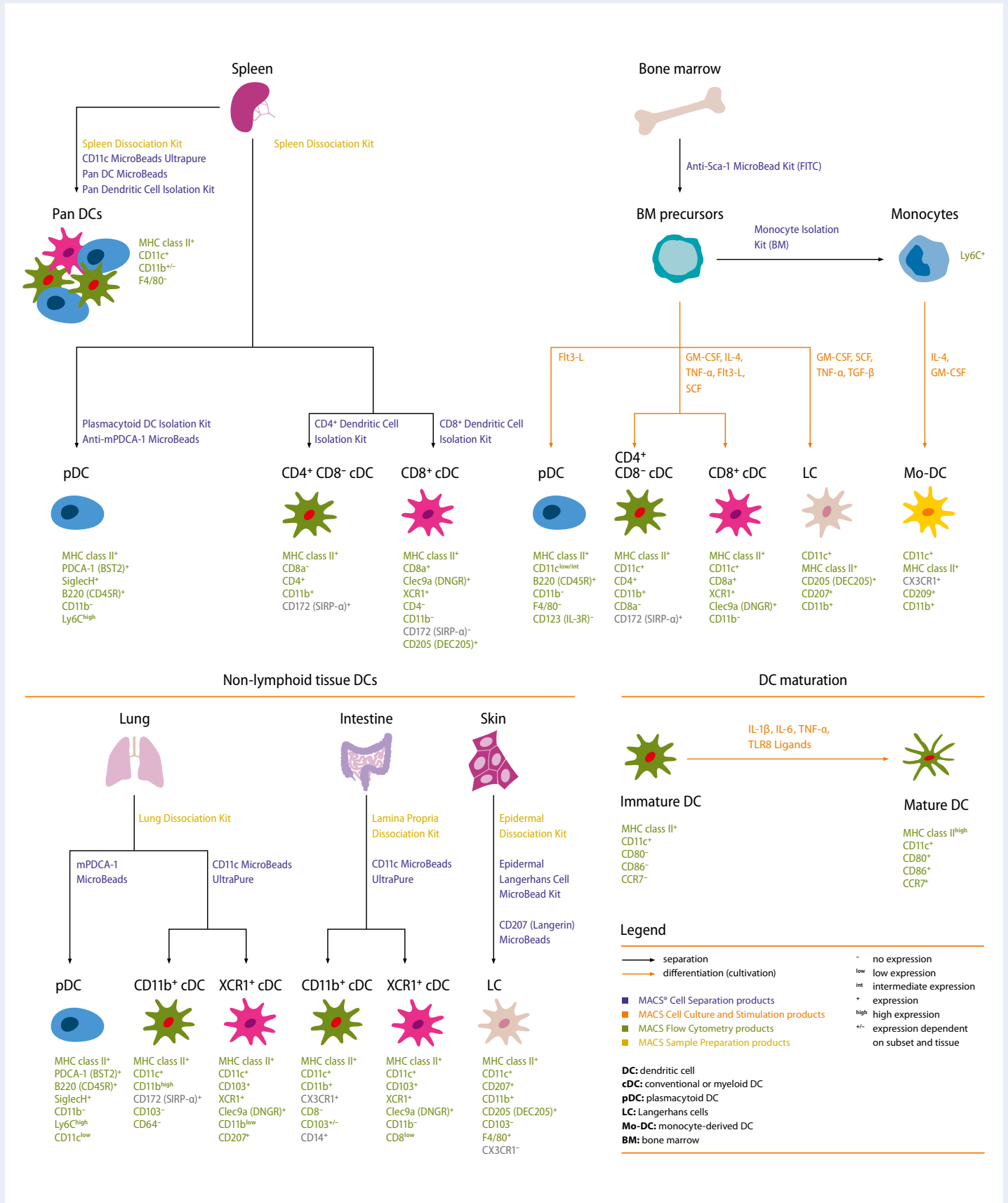
Products	Order no.
MC CD14 Monocyte Cocktail	130-092-859
Mo-DC Differentiation Inspector	130-093-567
Blood Dendritic Cell Enumeration Kit	130-091-086 130-091-324

### CliniMACS® Reagents

Products
CliniMACS® CD14 Reagent
CliniMACS CD304 (BDCA-4) Reagent
CliniMACS CD1c (BDCA-1)-Biotin

# Mouse dendritic cell generation

## Overview



# Mouse

## Product tables

### Tissue dissociation reagents

Product	Order no.
Spleen Dissociation Kit	130-095-926
Lung Dissociation Kit	130-095-927
Epidermis Dissociation Kit	130-095-928
Lamina Propria Dissociation Kit	130-097-410
Multi Tissue Dissociation Kit 1	130-110-201
Multi Tissue Dissociation Kit 2	130-110-203
Multi Tissue Dissociation Kit 3	130-110-204

For more information visit:

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

### Cell separation reagents

Product	Order no.
CD11c MicroBeads Ultrapure	130-108-338
Pan DC MicroBeads	130-092-465
Pan Dendritic Cell Isolation Kit	130-100-875
Plasmacytoid Dendritic Cell Isolation Kit	130-107-093
Anti-mPDCA-1 MicroBeads	130-091-965
CD4 <sup>+</sup> Dendritic Cell Isolation Kit	130-091-262
CD8 <sup>+</sup> Dendritic Cell Isolation Kit	130-091-169
Epidermal Langerhans Cell MicroBead Kit	130-095-408
CD207 (Langerin) MicroBeads	130-097-572

For more information visit:

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

### Cell culture reagents

Product	Research grade	Premium grade
Mouse IL-4	●	●
Mouse GM-CSF	●	●
Mouse Flt3-Ligand	●	●
Mouse TNF $\alpha$	●	●
Mouse IL-6	●	●
Mouse IL-1 $\beta$	●	●
Mouse SCF	●	●
Human TGF- $\beta$ 1		●
TLR9-Ligands	●	
TLR7/8-Ligands	●	

For more information about custom fillings and cocktails visit:

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



For more information about our instruments visit:

[www.miltenyibiotec.com/MACS-Automation](http://www.miltenyibiotec.com/MACS-Automation)

[www.miltenyibiotec.com/flow-cytometers](http://www.miltenyibiotec.com/flow-cytometers)

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

# Mouse

## Product tables

### Flow cytometry antibodies

Antigen	Vio-Blue®	Vio-Green®	FITC	VioBright™ FITC	PE	PerCP	PerCP-Vio® 700	PE-Vio 615	PE-Vio 770	APC	APC-Vio 770	Biotin	REAFinity
<b>Pan DC</b>													
MHC Class II	●	●	●		●		●			●		●	
CD11c	●		●		●		●		●	●	●	●	
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
F4/80			●		●		●	●	●	●		●	●
<b>pDC</b>													
MHC Class II	●	●	●		●		●			●		●	
CD11c	●		●		●		●		●	●	●	●	
mPDCA-1			●		●					●		●	
SiglecH			●		●					●		●	
CD45R (B220)	●	●	●		●	●	●		●	●	●	●	
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
Ly-6C	●	●	●		●		●		●	●	●	●	
<b>cDC</b>													
MHC Class II	●	●	●		●		●			●		●	
CD11c	●		●		●		●		●	●	●	●	
CD8a	●	●	●	●	●	●	●	●	●	●	●	●	●
CD4	●	●	●	●	●		●		●	●	●	●	●
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
CLEC9A				●	●							●	
XCR1				●	●	●	●		●	●	●	●	●
CD103			●		●					●		●	
CD64					●				●	●	●		●
CD205 (DEC205)					●				●	●		●	
CD207	●		●		●				●	●		●	
<b>LC</b>													
MHC Class II	●	●	●		●		●			●		●	
CD11c	●		●		●		●		●	●	●	●	
CD207	●		●		●				●	●		●	
CD11b	●	●	●	●	●		●	●	●	●	●	●	●
CD205 (DEC205)					●				●	●		●	
CD103			●		●					●		●	
F4/80			●		●		●	●	●	●	●	●	●

For more information about MACS Antibodies and recombinantly generated REAFinity Antibodies visit:

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

For customized filling or bulk requests of pure antibodies for mass cytometry please go to: [www.miltenyibiotec.com/CytoF](http://www.miltenyibiotec.com/CytoF)





## Miltenyi Biotec

### Germany/Austria/ Switzerland

Miltenyi Biotec GmbH  
Friedrich-Ebert-Straße 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 16A, 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.  
Schipholweg 68 H  
2316 XE 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 Biotech Technology &  
Trading (Shanghai) Co., Ltd.  
Room 2309  
No. 319, Xianxia Road  
Changning District  
200051 Shanghai, P.R. China  
Phone +86 21 62351005  
Fax +86 21 62350953  
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 Persicetana, 2/D  
40012 Calderara di Reno (BO)  
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

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

Miltenyi Biotec provides products and services worldwide. Visit [www.miltenyibiotec.com/local](http://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® GMP Products are for research use and ex vivo cell culture processing only, and are not intended for human in vivo applications. For regulatory status in the USA, please contact your local representative. MACS GMP Products are manufactured and tested under a quality system certified to ISO 13485 and are in compliance with relevant GMP guidelines. They are designed following the recommendations of USP <1043> on ancillary materials. The CliniMACS® System components, including Reagents, Tubing Sets, Instruments, and PBS/EDTA Buffer, are designed, manufactured and tested under a quality system certified to ISO 13485.

In the EU, the CliniMACS System components are available as CE-marked medical devices for their respective intended use, unless otherwise stated. The CliniMACS Reagents and Biotin Conjugates are intended for in vitro use only and are not designated for therapeutic use or direct infusion into patients. The CliniMACS Reagents in combination with the CliniMACS System are intended to separate human cells. Miltenyi Biotec as the manufacturer of the CliniMACS System does not give any recommendations regarding the use of separated cells for therapeutic purposes and does not make any claims regarding a clinical benefit. For the manufacturing and use of target cells in humans the national legislation and regulations - e.g. for the EU the Directive 2004/23/EC ("human tissues and cells"), or the Directive 2002/98/EC ("human blood and blood components") - must be followed. Thus, any clinical application of the target cells is exclusively within the responsibility of the user of a CliniMACS System.

In the US, the CliniMACS CD34 Reagent System, including the CliniMACS Plus Instrument, CliniMACS CD34 Reagent, CliniMACS Tubing Sets TS and LS, and the CliniMACS PBS/EDTA Buffer, is FDA approved; all other products of the CliniMACS Product Line are available for use only under an approved Investigational New Drug (IND) application or Investigational Device Exemption (IDE). CliniMACS MicroBeads are for research use only and not for human therapeutic or diagnostic use. CliniMACS, Vio, VioBright, VioBlue, VioGreen, MACS, and the MACS logo are registered trademarks or trademarks of Miltenyi Biotec GmbH and/or its affiliates in countries worldwide. Copyright © 2017 Miltenyi Biotec GmbH and/or its affiliates. All rights reserved.