

abcam

Immune cell markers

poster

References

References

1. Hoover-Plow J, G. Y. Challenges for heart disease stem cell therapy. *Vasc. Health Risk Manag.* 8, 99–113 (2012).
2. Warr, M. R., Pietras, E. M. & Passegué, E. Mechanisms controlling hematopoietic stem cell functions during normal hematopoiesis and hematological malignancies. *Wiley Interdiscip. Rev. Syst. Biol. Med.* 3, 681–701 (2011).
3. Weissman, I. L. & Shizuru, J. A. The origins of the identification and isolation of hematopoietic stem cells, and their capability to induce donor-specific transplantation tolerance and treat autoimmune diseases. *Blood* 112, 3543–3553 (2008).
4. Challen, G. a., Boles, N., Lin, K. K. & Goodell, M. a. Mouse Hematopoietic Stem Cell Identification and Analysis. *Cytom. A* 75, 14–24 (2009).
5. Sudo, T. et al. in *Stem Cell Biology in Normal Life and Diseases* (InTech, 2013). doi:10.5772/54474
6. Notta, F. et al. Isolation of Single Human Hematopoietic Stem Cells Capable of Long-Term Multilineage Engraftment. *Science* 333, 218–221 (2011).
7. Doulatov, S. et al. Revised map of the human progenitor hierarchy shows the origin of macrophages and dendritic cells in early lymphoid development. *Nat. Immunol.* 11, 585–93 (2010).
8. Majeti, R., Park, C. Y. & Weissman, I. L. Identification of a Hierarchy of Multipotent Hematopoietic Progenitors in Human Cord Blood. *Cell Stem Cell* 1, 635–645 (2007).
9. Goardon, N. et al. Coexistence of LMPP-like and GMP-like leukemia stem cells in acute myeloid leukemia. *Cancer Cell* 19, 138–152 (2011).
10. Welner, R. S., Pelayo, R. & Kincade, P. W. Evolving views on the genealogy of B cells. *Nat. Rev. Immunol.* 8, 95–106 (2008).
11. Doulatov, S., Notta, F., Laurenti, E. & Dick, J. E. Hematopoiesis: A human perspective. *Cell Stem Cell* 10, 120–136 (2012).
12. Mori, Y., Chen, J. Y., Pluvineau, J. V., Seita, J. & Weissman, I. L. Prospective isolation of human erythroid lineage-committed progenitors. *Proc. Natl. Acad. Sci.* 112, 9638–9643 (2015).
13. Chang, Y., Bluteau, D., Debili, N. & Vainchenker, W. From hematopoietic stem cells to platelets. *J. Thromb. Haemost.* 5, 318–327 (2007).
14. Novershtern, N. et al. Densely interconnected transcriptional circuits control cell states in human hematopoiesis. *Cell* 144, 296–309 (2011).
15. Lee, J. et al. Restricted dendritic cell and monocyte progenitors in human cord blood and bone marrow. *J. Exp. Med.* 212, 385–99 (2015).
16. Montaldo, E. et al. Human NK cell receptors/markers: A tool to analyze NK cell development, subsets and function. *Cytom. Part A* 83, 702–713 (2013).

17. Chen, Q. et al. Delineation of Natural Killer Cell Differentiation from Myeloid Progenitors in Human. *Sci. Rep.* 5, 15118 (2015).
18. Colucci, F., Caligiuri, M. a & Di Santo, J. P. What does it take to make a natural killer? *Nat. Rev. Immunol.* 3, 413–425 (2003).
19. Farag, S. S. & Caligiuri, M. A. Human natural killer cell development and biology. *Blood Rev.* 20, 123–137 (2006).
20. *Flow Cytometry.* (Humana Press, 2007). doi:10.1007/978-1-59745-451-3
21. Galon, J. et al. Type, density, and location of immune cells within human colorectal tumors predict clinical outcome. *Science* 313, 1960–4 (2006).
22. Finak, G. et al. Standardizing Flow Cytometry Immunophenotyping Analysis from the Human ImmunoPhenotyping Consortium. *Sci. Rep.* 6, 20686 (2016).
23. Kaminski, D. A., Wei, C., Qian, Y., Rosenberg, A. F. & Sanz, I. Advances in human B cell phenotypic profiling. *Front. Immunol.* 3, 1–15 (2012).
24. Orlic, D., Fischer, R., Nishikawa, S., Nienhuis, A. W. & Bodine, D. M. Purification and characterization of heterogeneous pluripotent hematopoietic stem cell populations expressing high levels of c-kit receptor. *Blood* 82, 762–70 (1993).
25. Bendall, S. C. et al. Single-Cell Trajectory Detection Uncovers Progression and Regulatory Coordination in Human B Cell Development. *Cell* 157, 714–725 (2014).
26. Wood, B. Multicolor Immunophenotyping: Human Immune System Hematopoiesis. *Methods Cell Biol.* 75, 559–576 (2004).
27. Yang, J., Zhang, L., Yu, C., Yang, X.-F. & Wang, H. Monocyte and macrophage differentiation: circulation inflammatory monocyte as biomarker for inflammatory diseases. *Biomark. Res.* 2, 1 (2014).
28. Murray, P. J. & Wynn, T. A. Protective and pathogenic functions of macrophage subsets. *Nat. Rev. Immunol.* 11, 723–37 (2011).
29. Pilling, D., Fan, T., Huang, D., Kaul, B. & Gomer, R. H. Identification of markers that distinguish monocyte-derived fibrocytes from monocytes, macrophages, and fibroblasts. *PLoS One* 4, 31–33 (2009).
30. Merad, M., Sathe, P., Helft, J., Miller, J. & Mortha, A. The dendritic cell lineage: ontogeny and function of dendritic cells and their subsets in the steady state and the inflamed setting. *Annu. Rev. Immunol.* 31, 563–604 (2013).
31. Elghetany, M. & Elghetany, M. Surface Antigen Changes during Normal Neutrophilic Development: A Critical Review. *Blood Cells, Mol. Dis.* 28, 260–274 (2002).
32. Mantovani, A., Cassatella, M. A., Costantini, C. & Jaillon, S. Neutrophils in the activation and regulation of innate and adaptive immunity. *Nat. Rev. Immunol.* 11, 519–531 (2011).

33. Behnen, M. et al. Immobilized Immune Complexes Induce Neutrophil Extracellular Trap Release by Human Neutrophil Granulocytes via Fc RIIIb and Mac-1. *J. Immunol.* 193, 1954–1965 (2014).
34. Lee, J. J. et al. Human versus mouse eosinophils: 'That which we call an eosinophil, by any other name would stain as red'. *J. Allergy Clin. Immunol.* 130, 572–584 (2012).
35. Mori, Y. et al. Identification of the human eosinophil lineage-committed progenitor: revision of phenotypic definition of the human common myeloid progenitor. *J. Exp. Med.* 206, 183–193 (2009).
36. Rothenberg, M. E. & Hogan, S. P. The eosinophil. *Annu. Rev. Immunol.* 24, 147–74 (2006).
37. Han, X. et al. Immunophenotypic study of basophils by multiparameter flow cytometry. *Arch. Pathol. Lab. Med.* 132, 813–819 (2008).
38. Cromheecke, J. L., Nguyen, K. T. & Huston, D. P. Emerging role of human basophil biology in health and disease. *Curr. Allergy Asthma Rep.* 14, 43–45 (2014).
39. Hauswirth, A. W. et al. Expression of cell surface antigens on mast cells: mast cell phenotyping. *Methods Mol. Biol.* 315, 77–90 (2006).
40. Sánchez-Muñoz, L., Teodósio, C., Morgado, J. M. & Escribano, L. Immunophenotypic characterization of bone marrow mast cells in mastocytosis and other mast cell disorders. *Methods Cell Biol.* 103, 333–59 (2011).
41. Murphy, K. *Janeway's Immunobiology.* (Garland Science, 2012).
42. Rodríguez-Perales, S. et al. Truncated RUNX1 protein generated by a novel t(1;21)(p32;q22) chromosomal translocation impairs the proliferation and differentiation of human hematopoietic progenitors. *Oncogene* 35, 1–10 (2015).
43. Deutsch, V. R. & Tomer, A. Megakaryocyte development and platelet production. *Br. J. Haematol.* 134, 453–466 (2006).
44. Van Velzen, J. F., Laros-Van Gorkom, B. A. P., Pop, G. A. M. & Van Heerde, W. L. Multicolor flow cytometry for evaluation of platelet surface antigens and activation markers. *Thromb. Res.* 130, 92–98 (2012).
45. Rector, K., Liu, Y. & Van Zant, G. Comprehensive hematopoietic stem cell isolation methods. *Methods Mol. Biol.* 976, 1–15 (2013).
46. Benveniste, P. et al. Intermediate-Term Hematopoietic Stem Cells with Extended but Time-Limited Reconstitution Potential. *Cell Stem Cell* 6, 48–58 (2010).
47. Qian, P. et al. The Dlk1-Gtl2 Locus Preserves LT-HSC Function by Inhibiting the PI3K-mTOR Pathway to Restrict Mitochondrial Metabolism. *Cell Stem Cell* 18, 214–28 (2015).
48. Oguro, H., Ding, L. & Morrison, S. J. SLAM family markers resolve functionally distinct subpopulations of hematopoietic stem cells and multipotent progenitors. *Cell Stem Cell* 13, 102–116 (2013).

49. Lebert-Ghali, C. É. et al. Hoxa cluster genes determine the proliferative activity of adult mouse hematopoietic stem and progenitor cells. *Blood* 127, 87–90 (2016).
50. Ema, H., Morita, Y. & Suda, T. Heterogeneity and hierarchy of hematopoietic stem cells. *Exp. Hematol.* 42, 74–82 (2014).
51. Ichim, C. V, Atkins, H. L., Iscove, N. N. & Wells, R. a. Identification of a role for the nuclear receptor EAR-2 in the maintenance of clonogenic status within the leukemia cell hierarchy. *Leuk. Off. J. Leuk. Soc. Am. Leuk. Res. Fund, U.K* 25, 1687–96 (2011).
52. Mantel, C. R. et al. Enhancing Hematopoietic Stem Cell Transplantation Efficacy by Mitigating Oxygen Shock. *Cell* 161, 1553–1565 (2015).
53. Chen, E. et al. Distinct effects of concomitant Jak2V617F expression and Tet2 loss in mice promote disease progression in myeloproliferative neoplasms. *Blood* 125, 327–335 (2015).
54. Cheng, C. W. et al. Prolonged fasting reduces IGF-1/PKA to promote hematopoietic-stem-cell- based regeneration and reverse immunosuppression. *Cell Stem Cell* 14, 810–823 (2014).
55. Takubo, K. et al. Regulation of glycolysis by Pdk functions as a metabolic checkpoint for cell cycle quiescence in hematopoietic stem cells. *Cell Stem Cell* 12, 49–61 (2013).
56. Miyawaki, K. et al. CD41 Marks the Initial Myelo-Erythroid Lineage Specification in Adult Mouse Hematopoiesis: Redefinition of Murine Common Myeloid Progenitor. *Stem Cells* 33, 976–987 (2015).
57. Yamamoto, R. et al. Clonal analysis unveils self-renewing lineage-restricted progenitors generated directly from hematopoietic stem cells. *Cell* 154, 1112–1126 (2013).
58. Yu, J., Freud, A. G. & Caligiuri, M. A. Location and cellular stages of natural killer cell development. *Trends Immunol.* 34, 573–582 (2013).
59. Klose, C. S. N. et al. Differentiation of type 1 ILCs from a common progenitor to all helper-like innate lymphoid cell lineages. *Cell* 157, 340–356 (2014).
60. Nakorn, T. N., Miyamoto, T. & Weissman, I. L. Characterization of mouse clonogenic megakaryocyte progenitors. *Proc. Natl. Acad. Sci. U. S. A.* 100, 205–210 (2003).
61. Huntington, N. D., Vosshenrich, C. a J. & Di Santo, J. P. Developmental pathways that generate natural-killer-cell diversity in mice and humans. *Nat. Rev. Immunol.* 7, 703–14 (2007).
62. Yui, M. A. & Rothenberg, E. V. Developmental gene networks: a triathlon on the course to T cell identity. *Nat. Rev. Immunol.* 14, 529–45 (2014).
63. Mukai, K. et al. Critical role of P1-Runx1 in mouse basophil development. *Blood* 120, 76–85 (2012).
64. Terry, R. L. & Miller, S. D. Molecular control of monocyte development. *Cell. Immunol.* 291, 16–21 (2014).

65. Tamoutounour, S. et al. Origins and Functional Specialization of Macrophages and of Conventional and Monocyte-Derived Dendritic Cells in Mouse Skin. *Immunity* 39, 925–938 (2013).
66. Misharin, A. V., Morales-Nebreda, L., Mutlu, G. M., Budinger, G. R. S. & Perlman, H. Flow cytometric analysis of macrophages and dendritic cell subsets in the mouse lung. *Am. J. Respir. Cell Mol. Biol.* 49, 503–510 (2013).
67. Gautier, E. L. et al. Gene-expression profiles and transcriptional regulatory pathways that underlie the identity and diversity of mouse tissue macrophages. *Nat. Immunol.* 13, 1118–28 (2012).
68. Imai, Y. et al. Skin-specific expression of IL-33 activates group 2 innate lymphoid cells and elicits atopic dermatitis-like inflammation in mice. *Proc. Natl. Acad. Sci. U. S. A.* 110, 13921–6 (2013).
69. Christoffersson, G. et al. VEGF-A recruits a proangiogenic MMP-9-delivering neutrophil subset that induces angiogenesis in transplanted hypoxic tissue. *Blood* 120, 4653–4662 (2012).
70. Cho, H. et al. The Loss of RGS Protein-G i2 Interactions Results in Markedly Impaired Mouse Neutrophil Trafficking to Inflammatory Sites. *Mol. Cell. Biol.* 32, 4561–4571 (2012).
71. Rose, S., Misharin, A. & Perlman, H. A novel Ly6C/Ly6G-based strategy to analyze the mouse splenic myeloid compartment. *Cytom. Part A* 81 A, 343–350 (2012).
72. Doyle, A. D. et al. Expression of the secondary granule proteins major basic protein 1 (MBP-1) and eosinophil peroxidase (EPX) is required for eosinophilopoiesis in mice. *Blood* 122, 781–790 (2013).
73. Iwamoto, H., Matsubara, T., Nakazato, Y., Namba, K. & Takeda, Y. Decreased expression of CD200R3 on mouse basophils as a novel marker for IgG1-mediated anaphylaxis. *Immunity, Inflamm. Dis.* 3, 280–288 (2015).
74. Obata, K. et al. Basophils are essential initiators of a novel type of chronic allergic inflammation. *Blood* 110, 913–920 (2007).
75. Bakocevic, N. et al. CD41 is a reliable identification and activation marker for murine basophils in the steady state and during helminth and malarial infections. *Eur. J. Immunol.* 44, 1823–1834 (2014).
76. Dahlin, J. S. & Hallgren, J. Mast cell progenitors: Origin, development and migration to tissues. *Mol. Immunol.* 63, 9–17 (2015).
77. S., D., DingZhoujie & HallgrenJenny. Distinguishing Mast Cell Progenitors from Mature Mast Cells in Mice. *Stem Cells Dev.* 24, 1703–1711 (2015).
78. Chen, C. C., Grimbaldston, M. A., Tsai, M., Weissman, I. L. & Galli, S. J. Identification of mast cell progenitors in adult mice. *Proc. Natl. Acad. Sci. U. S. A.* 102, 11408–13 (2005).
79. Gri, G. et al. Mast cell: An emerging partner in immune interaction. *Front. Immunol.* 3, 1–11 (2012).

80. Garcia-Santos, D. et al. Heme oxygenase 1 is expressed in murine erythroid cells where it controls the level of regulatory heme. *Blood* 123, 2269–2277 (2014).
81. Cuyper, I. M. De et al. A novel flow cytometry – based platelet aggregation assay. *Blood* 121, 70–80 (2013).
82. Woolthuis, C. M. & Park, C. Y. Hematopoietic stem/progenitor cell commitment to the megakaryocyte lineage. *Blood* 127, 1242–1249 (2016).
83. Niswander, L. M., McGrath, K. E., Kennedy, J. C. & Palis, J. Improved quantitative analysis of primary bone marrow megakaryocytes utilizing imaging flow cytometry. *Cytom. Part A* 85, 302–312 (2014).
84. Pronk, C. J. H. et al. Elucidation of the Phenotypic, Functional, and Molecular Topography of a Myeloerythroid Progenitor Cell Hierarchy. *Cell Stem Cell* 1, 428–442 (2007).
85. Yu, M. & Cantor, A. B. Megakaryopoiesis and thrombopoiesis: an update on cytokines and lineage surface markers. *Methods Mol. Biol.* 788, 291–303 (2012).
86. Boilard, E. et al. Platelets Amplify Inflammation in Arthritis via Collagen-Dependent Microparticle Production. *Science* 327, 580–583 (2010).
87. Moser, M., Nieswandt, B., Ussar, S., Pozgajova, M. & Fässler, R. Kindlin-3 is essential for integrin activation and platelet aggregation. *Nat. Med.* 14, 325–330 (2008).
88. Nieswandt, B. et al. Loss of talin1 in platelets abrogates integrin activation, platelet aggregation, and thrombus formation in vitro and in vivo. *J. Exp. Med.* 204, 3113–8 (2007).
89. Cabezas-Wallscheid, N. & Trumpp, A. Potency finds its niches. *Science INSIGHTS | Perspectives* 351, 126–127 (2016).
90. Notta, F. et al. Distinct routes of lineage development reshape the human blood hierarchy across ontogeny. *Science* 351, 1–16 (2015).
91. Paul, F. et al. Transcriptional Heterogeneity and Lineage Commitment in Myeloid Progenitors. *Cell* 163, 1663–1677 (2015).

Other research articles reviewed

- Abeles, R. D. et al. CD14, CD16 and HLA-DR reliably identifies human monocytes and their subsets in the context of pathologically reduced HLA-DR expression by CD14hi/CD16neg monocytes: Expansion of CD14hi/CD16pos and contraction of CD14lo/CD16pos monocytes in acute liver fail. *Cytom. Part A* 81 A, 823–834 (2012).
- Adolfsson, J. et al. Identification of Flt3+ lympho-myeloid stem cells lacking erythromegakaryocytic potential: A revised road map for adult blood lineage commitment. *Cell* 121, 295–306 (2005).
- Ambarus, C. A. et al. Systematic validation of specific phenotypic markers for in vitro polarized human macrophages. *J. Immunol. Methods* 375, 196–206 (2012).
- Amsen, D., Helbig, C. & Backer, R. A. Notch in T Cell Differentiation: All Things Considered. *Trends Immunol.* 36, 802–814 (2015).
- Anderson, S. M., Tomayko, M. M., Ahuja, A., Haberman, A. M. & Shlomchik, M. J. New markers for murine memory B cells that define mutated and unmutated subsets. *J. Exp. Med.* 204, 2103–14 (2007).
- Anjos-Afonso, F. et al. CD34-cells at the apex of the human hematopoietic stem cell hierarchy have distinctive cellular and molecular signatures. *Cell Stem Cell* 13, 161–174 (2013).
- Ardavín, C. Origin, precursors and differentiation of mouse dendritic cells. *Nat. Rev. Immunol.* 3, 582–590 (2003).
- Auffray, C. et al. CX3CR1+ CD115+ CD135+ common macrophage/DC precursors and the role of CX3CR1 in their response to inflammation. *J. Exp. Med.* 206, 595–606 (2009).
- Barteneva, N. S., Editors, I. A. V. & Walker, J. M. *Imaging Flow Cytometry*. (Springer New York, 2016). doi:10.1007/978-1-4939-3302-0
- Baumgarth, N. & Roederer, M. A practical approach to multicolor flow cytometry for immunophenotyping. *J. Immunol. Methods* 243, 77–97 (2000).
- Bell, K. S. et al. The role of individual protein kinase C isoforms in mouse mast cell function and their targeting by the immunomodulatory parasitic worm product, ES-62. *Immunol. Lett.* 168, 31–40 (2015).
- Belz, G. T. & Nutt, S. L. Transcriptional programming of the dendritic cell network. *Nat. Rev. Immunol.* 12, 101–13 (2012).
- Bendall, S. C. et al. Single-Cell Mass Cytometry of Differential Immune and Drug Responses Across a Human Hematopoietic Continuum. *Science* 332, 687–696 (2011).
- Bianchi, E., Norfo, R., Pennucci, V., Zini, R. & Manfredini, R. Genomic landscape of megakaryopoiesis and platelet function defects. *Blood* 127, 1249–1260 (2016).

- Blasi, T. et al. Label-free cell cycle analysis for high-throughput imaging flow cytometry. *Nat. Commun.* 7, 10256 (2016).
- Bleesing, J. J. H. & Fleisher, T. a. Human B cells express a CD45 isoform that is similar to murine B220 and is downregulated with acquisition of the memory B-cell marker CD27. *Cytometry. Part B, Clinical cytometry* 51, 1–8 (2003).
- Blom, B. & Spits, H. Development of Human Lymphoid Cells. *Annu. Rev. Immunol.* 24, 287–320 (2006).
- T-Cell Development. (Springer New York, 2016). doi: 10.1007/978-1-4939-2809-5
- Boyer, S. W., Schroeder, A. V., Smith-Berdan, S. & Forsberg, E. C. All Hematopoietic Cells Develop from Hematopoietic Stem Cells through Flk2/Flt3-Positive Progenitor Cells. *Cell Stem Cell* 9, 64–73 (2011).
- Cabezas-Wallscheid, N. & Trumpp, A. STEM CELLS. Potency finds its niches. *Science* 351, 126-7. (2016).
- Calderón-Gómez, E. & Fillatreau, S. Regulatory B Cells. *Methods Mol. Biol.* 1190, 105–13 (2014).
- Cassado, A. dos A., D'Império Lima, M. R. & Bortoluci, K. R. Revisiting Mouse Peritoneal Macrophages: Heterogeneity, Development, and Function. *Front. Immunol.* 6, 225 (2015).
- Cassard, L., Jonsson, F., Arnaud, S. & Daeron, M. Fc Receptors Inhibit Mouse and Human Basophil Activation. *J. Immunol.* 189, 2995–3006 (2012).
- Ceredig, R., Rolink, A. G. & Brown, G. Models of haematopoiesis: seeing the wood for the trees. *Nat. Rev. Immunol.* 9, 293–300 (2009).
- Chen, Z., Hu, M. & Shivdasani, R. A. Expression analysis of primary mouse megakaryocyte differentiation and its application in identifying stage-specific molecular markers and a novel transcriptional target of NF-E2. *Blood* 109, 1451–9 (2007).
- Chotinantakul, K. & Leraanansaksiri, W. Hematopoietic Stem Cell Development, Niches, and Signaling Pathways. *Bone Marrow Res.* 2012, 1–16 (2012).
- Chuluundorj, D., Harding, S. A., Abernethy, D. & La Flamme, A. C. Expansion and preferential activation of the CD14(+)CD16(+) monocyte subset during multiple sclerosis. *Immunol. Cell Biol.* 92, 509–517 (2014).
- Collison, J. L., Carlin, L. M., Eichmann, M., Geissmann, F. & Peakman, M. Heterogeneity in the Locomotory Behavior of Human Monocyte Subsets over Human Vascular Endothelium In Vitro. *J. Immunol.* 195, 1162–1170 (2015).
- Conley, J. M., Gallagher, M. P. & Berg, L. J. T Cells and Gene Regulation: The Switching On and Turning Up of Genes after T Cell Receptor Stimulation in CD8 T Cells. *Front. Immunol.* 7, 1–7 (2016).
- Dalli, J. et al. The novel 13S,14S-epoxy-maresin converted by human macrophages to maresin 1 (MaR1), inhibits leukotriene A4 hydrolase (LTA4H), and shifts macrophage phenotype. *FASEB J.* 27, 2573–2583 (2013).

- Davies, L. C., Jenkins, S. J., Allen, J. E. & Taylor, P. R. Tissue-resident macrophages. *Nat. Immunol.* 14, 986–995 (2013).
- Dawicki, W. & Marshall, J. S. New and emerging roles for mast cells in host defence. *Curr. Opin. Immunol.* 19, 31–38 (2007).
- De Rosa, S. C., Herzenberg, L. a, Herzenberg, L. a & Roederer, M. 11-color, 13-parameter flow cytometry: identification of human naive T cells by phenotype, function, and T-cell receptor diversity. *Nat. Med.* 7, 245–248 (2001).
- De Rosa, S. C., Brenchley, J. M. & Roederer, M. Beyond six colors: a new era in flow cytometry. *Nat. Med.* 9, 112–117 (2003).
- Dong, Z. M. & Murphy, J. W. Cryptococcal polysaccharides induce L-selectin shedding and tumor necrosis factor receptor loss from the surface of human neutrophils. *J. Clin. Invest.* 97, 689–698 (1996).
- Duan, M. et al. CD11b immunophenotyping identifies inflammatory profiles in the mouse and human lungs. *Mucosal Immunol.* 9, 1–14 (2015).
- DuMont, A. L. et al. Staphylococcus aureus LukAB cytotoxin kills human neutrophils by targeting the CD11b subunit of the integrin Mac-1. *Proc. Natl. Acad. Sci. U. S. A.* 110, 10794–9 (2013).
- Durafourt, B. A. et al. Comparison of polarization properties of human adult microglia and blood-derived macrophages. *Glia* 60, 717–727 (2012).
- Dvorak, A. M. The mouse basophil, a rare and rarely recognized granulocyte. *Blood* 96, 1616–7 (2000).
- Ellington, E. H. et al. Using multiple imputation to estimate missing data in meta-regression. *Methods in Ecology and Evolution* 6, (2015).
- Ellmeier, W., Haust, L. & Tschismarov, R. Transcriptional control of CD4 and CD8 coreceptor expression during T cell development. *Cell. Mol. Life Sci.* 70, 4537–4553 (2013).
- Erwig, L. P. & Gow, N. A. R. Interactions of fungal pathogens with phagocytes. *Nat. Rev. Microbiol.* 14, 163–176 (2016).
- Ethier, C., Lacy, P. & Davoine, F. Identification of human eosinophils in whole blood by flow cytometry. *Methods Mol. Biol.* 1178, 81–92 (2014).
- Falcone, F. H., Haas, H. & Gibbs, B. F. The human basophil: a new appreciation of its role in immune responses. *Blood* 96, 4028–38 (2000).
- Fares, I. et al. Pyrimidoindole derivatives are agonists of human hematopoietic stem cell self-renewal. *Science* 345, 1509 (2014).
- Ferreirós-Vidal, I. et al. Genome-wide identification of Ikaros targets elucidates its contribution to mouse B-cell lineage specification and pre-B-cell differentiation. *Blood* 121, 1769–82 (2013).

- Franco, C. B., Chen, C. C., Drukker, M., Weissman, I. L. & Galli, S. J. Distinguishing Mast Cell and Granulocyte Differentiation at the Single-Cell Level. *Cell Stem Cell* 6, 361–368 (2010).
- Freitas, M., Porto, G., Lima, J. L. F. C. & Fernandes, E. Isolation and activation of human neutrophils in vitro. The importance of the anticoagulant used during blood collection. *Clin. Biochem.* 41, 570–575 (2008).
- Frelin, C. et al. GATA-3 regulates the self-renewal of long-term hematopoietic stem cells. *Nat. Immunol.* 14, 1037–44 (2013).
- Freud, A. G. & Caligiuri, M. A. Human natural killer cell development. *Immunol. Rev.* 214, 56–72 (2006).
- Freud, A. G., Yu, J. & Caligiuri, M. A. Human natural killer cell development in secondary lymphoid tissues. *Semin. Immunol.* 26, 132–137 (2014).
- Fu, B., Tian, Z. & Wei, H. Subsets of human natural killer cells and their regulatory effects. *Immunology* 141, 483–489 (2014).
- Fulkerson, P. C. & Rothenberg, M. E. Targeting eosinophils in allergy, inflammation and beyond. *Nat. Rev. Drug Discov.* 12, 117–29 (2013).
- Gay, L. J. & Felding-Habermann, B. Contribution of platelets to tumour metastasis. *Nat. Rev. Cancer* 11, 123–34 (2011).
- Geissman, F. et al. Development of monocytes, macrophages, and dendritic cells. *J. Microbiol. Immunol. Infect.* 656–662 (2010).
- Germain, R. N. T-cell development and the CD4-CD8 lineage decision. *Nat. Rev. Immunol.* 2, 309–322 (2002).
- Giepmans, B., Adams, S., Ellisman, M. & Tsien, R. The Fluorescent Toolbox for Assessing Protein Location and Function. *Science* 217–224 (2006).
- Gold, M., Marsolais, D. & Blanchet, M.-R. Mast Cells - Methods and Protocols. *Methods in Molecular Biology* 1220, (2015).
- Gordon, S., Plüddemann, A. & Martinez Estrada, F. Macrophage heterogeneity in tissues: Phenotypic diversity and functions. *Immunol. Rev.* 262, 36–55 (2014).
- Gordon, S. & Taylor, P. R. Monocyte and macrophage heterogeneity. *Nat. Rev. Immunol.* 5, 953–964 (2005).
- Guilliams, M. et al. Dendritic cells, monocytes and macrophages: a unified nomenclature based on ontogeny. *Nat. Rev. Immunol.* 14, 571–8 (2014).
- Gurish, M. F. & Austen, K. F. Developmental Origin and Functional Specialization of Mast Cell Subsets. *Immunity* 37, 25–33 (2012).
- Haley, K. M., Recht, M. & McCarty, O. J. T. Neonatal platelets: mediators of primary hemostasis in the developing hemostatic system. *Pediatr. Res.* 76, 230–7 (2014).

- Hayakawa, Y., Huntington, N. D., Nutt, S. L. & Smyth, M. J. Functional subsets of mouse natural killer cells. *Immunol. Rev.* 214, 47–55 (2006).
- Heazlewood, S. Y., Williams, B., Storan, M. J. & Nilsson, S. K. The prospective isolation of viable, high ploidy megakaryocytes from adult murine bone marrow by fluorescence activated cell sorting. *Methods Mol. Biol.* 1035, 121–133 (2013).
- Heib, V. et al. Mast cells are crucial for early inflammation, migration of Langerhans cells, and CTL responses following topical application of TLR7 ligand in mice. *Blood* 110, 946–953 (2007).
- Hendrix, S. et al. Mast cells protect from post-traumatic brain inflammation by the mast cell-specific chymase mouse mast cell protease-4. *FASEB J.* 27, 920–929 (2013).
- Herzenberg, L. A., Tung, J., Moore, W. A., Herzenberg, L. A. & Parks, D. R. Interpreting flow cytometry data: a guide for the perplexed. *Nat. Immunol.* 7, 681–5 (2006).
- Herzog, S., Reth, M. & Jumaa, H. Regulation of B-cell proliferation and differentiation by pre-B-cell receptor signalling. *Nat. Rev. Immunol.* 9, 195–205 (2009).
- Hess, D. C. et al. Hematopoietic origin of microglial and perivascular cells in brain. *Exp. Neurol.* 186, 134–144 (2004).
- Hirahara, K. & Nakayama, T. CD4 + T-cell subsets in inflammatory diseases: beyond the Th1/Th2 paradigm. *Int. Immunol.* 28, 163–171 (2016).
- Holm, M. et al. Seven week culture of functional human mast cells from buffy coat preparations. *J. Immunol. Methods* 336, 213–221 (2008).
- Huang, H. & Li, Y. Mechanisms controlling mast cell and basophil lineage decisions. *Curr. Allergy Asthma Rep.* 14, 457 (2014).
- Hulspas, R., O’Gorman, M. R. G., Wood, B. L., Gratama, J. W. & Robert Sutherland, D. Considerations for the control of background fluorescence in clinical flow cytometry. *Cytom. Part B - Clin. Cytom.* 76, 355–364 (2009).
- Ingersoll, M. et al. Comparison of gene expression profiles between human and mouse monocyte subsets. *Blood* 115, 10–20 (2010).
- Ivanova, N. B. et al. A stem cell molecular signature. *Science* 298, 601–4 (2002).
- Jablonski, K. A. et al. Novel Markers to Delineate Murine M1 and M2 Macrophages. *PLoS One* 10, e0145342 (2015).
- Jensen, C. T., Strid, T. & Sigvardsson, M. Exploring the multifaceted nature of the common lymphoid progenitor compartment. *Curr. Opin. Immunol.* 39, 121–126 (2016).
- Josefowicz, S. Z., Lu, L. F. & Rudensky, A. Y. Regulatory T cells: mechanisms of differentiation and function. *Annu. Rev. Immunol.* 30, 531–564 (2012).
- Kawamura, S., Onai, N., Takenaka, K., Akashi, K. & Ohteki, T. Identification of common monocyte progenitors, pre-monocytes, and granulocyte monocyte progenitors in human umbilical cord blood. *Exp. Hematol.* 43, S72 (2015).

- Khan, J. A. et al. Fetal liver hematopoietic stem cell niches associate with portal vessels. *Science* 351, 176–180 (2016).
- Kiel, M. J. et al. SLAM family receptors distinguish hematopoietic stem and progenitor cells and reveal endothelial niches for stem cells. *Cell* 121, 1109–1121 (2005).
- Kigerl, K. A. et al. Identification of Two Distinct Macrophage Subsets with Divergent Effects Causing either Neurotoxicity or Regeneration in the Injured Mouse Spinal Cord. *J. Neurosci.* 29, 13435–13444 (2009).
- Kina, T. et al. The monoclonal antibody TER-119 recognizes a molecule associated with glycophorin A and specifically marks the late stages of murine erythroid lineage. *Br. J. Haematol.* 109, 280–287 (2000).
- Kirshenbaum, A. S. et al. Demonstration that human mast cells arise from a progenitor cell population that is CD34+, c-kit+, and expresses aminopeptidase N (CD13). *Blood* 94, 2333–2342 (1999).
- Klimchenko, O. et al. A common bipotent progenitor generates the erythroid and megakaryocyte lineages in embryonic stem cell-derived primitive hematopoiesis. *Blood* 114, 1506–1517 (2009).
- Knol, E. F. & Gibbs, B. F. Basophils and antigen presentation: Of mice and not men? *Allergy Eur. J. Allergy Clin. Immunol.* 67, 579–580 (2012).
- Kobayashi, H., Ogawa, M., Alford, R., Choyke, P. L. & Urano, Y. New Strategies for Fluorescent Probe Design in Medical Diagnostic Imaging. *Chem. Rev.* 110, 2620–2640 (2010).
- Koch, U. & Radtke, F. Mechanisms of T Cell Development and Transformation. *Annu. Rev. Cell Dev. Biol.* 27, 539–562 (2011).
- Kochenderfer, J. N. & Rosenberg, S. A. Treating B-cell cancer with T cells expressing anti-CD19 chimeric antigen receptors. *Nat. Rev. Clin. Oncol.* 10, 267–276 (2013).
- Kolaczowska, E. & Kubes, P. Neutrophil recruitment and function in health and inflammation. *Nat. Rev. Immunol.* 13, 159–75 (2013).
- Koyanagi, A., Sekine, C. & Yagita, H. Expression of Notch receptors and ligands on immature and mature T cells. *Biochem. Biophys. Res. Commun.* 418, 799–805 (2012).
- Mast Cells. (Humana Press, 2015). doi: 10.1007/978-1-4939-1568-2.
- Kurosaki, T., Shinohara, H. & Baba, Y. B cell signaling and fate decision. *Annu. Rev. Immunol.* 28, 21–55 (2010).
- Lai, L., Alaverdi, N., Maltais, L. & Morse, H. C. Mouse cell surface antigens: nomenclature and immunophenotyping. *J. Immunol.* 160, 3861–8 (1998).
- Lämmermann, T. et al. Neutrophil swarms require LTB4 and integrins at sites of cell death in vivo. *Nature* 498, 371–5 (2013).
- Lavin, Y., Mortha, A., Rahman, A. & Merad, M. Regulation of macrophage development and function in peripheral tissues. *Nat. Rev. Immunol.* 15, 731–44 (2015).

- Lee, J. J. & McGarry, M. P. When is a mouse basophil not a basophil? *Blood* 109, 859–861 (2006).
- Loh, Y. H. et al. Reprogramming of T cells from human peripheral blood. *Cell Stem Cell* 7, 15–19 (2010).
- Maecker, H. T. & Trotter, J. Flow cytometry controls, instrument setup, and the determination of positivity. *Cytom. Part A* 69A, 1037–1042 (2006).
- Mantovani, A. et al. The chemokine system in diverse forms of macrophage activation and polarization. *Trends Immunol.* 25, 677–686 (2004).
- Martinez, F. O. & Gordon, S. The M1 and M2 paradigm of macrophage activation: time for reassessment. *F1000Prime Rep.* 6, 13 (2014).
- Martinez, F. O., Gordon, S., Locati, M. & Mantovani, A. Transcriptional Profiling of the Human Monocyte-to-Macrophage Differentiation and Polarization: New Molecules and Patterns of Gene Expression. *J. Immunol.* 177, 7303–7311 (2006).
- Matsumura-Takeda, K. et al. CD41+/CD45+ cells without acetylcholinesterase activity are immature and a major megakaryocytic population in murine bone marrow. *Stem Cells* 25, 862–70 (2007).
- Metcalfe, D. D. Mast cells and mastocytosis. *Blood* 112, 946–956 (2008).
- Mommert, S. et al. Human basophil chemotaxis and activation are regulated via the histamine H4 receptor. [Published online ahead of print March 10 2016]. *Allergy* (2016). doi:10.1111/all.12875
- Morrison, S. J., Wandycz, A. M., Akashi, K., Globerson, A. & Weissman, I. L. The aging of hematopoietic stem cells. *Nat. Med.* 2, 1011–6 (1996).
- Morrison, S. J. & Weissman, I. L. The long-term repopulating subset of hematopoietic stem cells is deterministic and isolatable by phenotype. *Immunity* 1, 661–673 (1994).
- Munoz, N. M. & Leff, A. R. Highly purified selective isolation of eosinophils from human peripheral blood by negative immunomagnetic selection. *Nat. Protoc.* 1, 2613–2620 (2006).
- Murray, P. J. et al. Macrophage Activation and Polarization: Nomenclature and Experimental Guidelines. *Immunity* 41, 14–20 (2014).
- Nacionales, D. C. et al. Aged Mice Are Unable to Mount an Effective Myeloid Response to Sepsis. *J. Immunol.* 192, 612–622 (2014).
- Nelissen, S. et al. Mast cells protect from post-traumatic spinal cord damage in mice by degrading inflammation-associated cytokines via mouse mast cell protease 4. *Neurobiol. Dis.* 62, 260–272 (2014).
- Nguyen, J. et al. Morphine stimulates cancer progression and mast cell activation and impairs survival in transgenic mice with breast cancer. *Br. J. Anaesth.* 113, 1–10 (2014).

- Notta, F. et al. Distinct routes of lineage development reshape the human blood hierarchy across ontogeny. *Science* 351, 1–16 (2015).
- O'Donnell, E. a, Ernst, D. N. & Hingorani, R. Multiparameter flow cytometry: advances in high resolution analysis. *Immune Netw.* 13, 43–54 (2013).
- Parekh, C. & Crooks, G. M. Critical differences in hematopoiesis and lymphoid development between humans and mice. *J. Clin. Immunol.* 33, 711–715 (2013).
- Paul, F. et al. Transcriptional Heterogeneity and Lineage Commitment in Myeloid Progenitors. *Cell* 163, 1663–1677 (2015).
- Perdiguerro, E. G. & Geissmann, F. The development and maintenance of resident macrophages. *Nat. Immunol.* 17, 2–8 (2015).
- Perfetto, S. P. et al. Amine reactive dyes: An effective tool to discriminate live and dead cells in polychromatic flow cytometry. *J. Immunol. Methods* 313, 199–208 (2006).
- Perfetto, S. P., Chattopadhyay, P. K. & Roederer, M. Seventeen-colour flow cytometry: unravelling the immune system. *Nat. Rev. Immunol.* 4, 648–55 (2004).
- Piliponsky, A. M. et al. The chymase mouse mast cell protease 4 degrades TNF, limits inflammation, and promotes survival in a model of sepsis. *Am. J. Pathol.* 181, 875–886 (2012).
- Punt, J., Owen, J. & Caligiuri, M. A. The biology of human natural killer-cell subsets. *Trends Immunol.* 22, 633–640 (2001).
- Radtke, F., MacDonald, H. R. & Tacchini-Cottier, F. Regulation of innate and adaptive immunity by Notch. *Nat. Rev. Immunol.* 13, 427–37 (2013).
- Ribatti, D. & Crivellato, E. in *Mast Cells and Tumours* 3–48 (Springer Netherlands, 2011). doi:10.1007/978-94-007-1469-4_2
- Roederer, M. Spectral compensation for flow cytometry: visualization artifacts, limitations, and caveats. *Cytometry* 45, 194–205 (2001).
- Rogers, N. M., Ferenbach, D. a, Isenberg, J. S., Thomson, A. W. & Hughes, J. Dendritic cells and macrophages in the kidney: a spectrum of good and evil. *Nat. Rev. Nephrol.* 10, 625–43 (2014).
- Rosser, E. C. & Mauri, C. Regulatory B cells in experimental mouse models of arthritis. *Methods in molecular biology* (Clifton, N.J.) 1190, (2014).
- Rothenberg, E. V. & Taghon, T. Molecular Genetics of T Cell Development. *Annu. Rev. Immunol.* 23, 601–649 (2005).
- Rousseau, M. et al. Detection and quantification of microparticles from different cellular lineages using flow cytometry. Evaluation of the impact of secreted phospholipase A2 on microparticle assessment. *PLoS One* 10, 1–27 (2015).
- Schulze, H. Culture, Expansion, and Differentiation of Murine Megakaryocytes from Fetal Liver, Bone Marrow, and Spleen. *Curr. Protoc. Immunol.* 22F.6.1–22F.6.15 (2016).

- Seki, T. et al. Generation of induced pluripotent stem cells from human terminally differentiated circulating t cells. *Cell Stem Cell* 7, 11–13 (2010).
- Sellge, G. & Bischoff, S. C. Isolation, culture, and characterization of intestinal mast cells. *Methods Mol. Biol.* 315, 123–138 (2006).
- Semple, J. W., Italiano, J. E. & Freedman, J. Platelets and the immune continuum. *Nat. Rev. Immunol.* 11, 264–74 (2011).
- Seo, W. & Taniuchi, I. Transcriptional regulation of early T-cell development in the thymus. *Eur. J. Immunol.* 46, 531–538 (2016).
- Shi, G.-P., Bot, I. & Kovanen, P. T. Mast cells in human and experimental cardiometabolic diseases. *Nat. Rev. Cardiol.* 12, 643–58 (2015).
- Shortman, K. & Liu, Y.-J. Mouse and Human Dendritic Cell Subtypes. *Nat. Rev. Immunol.* 2, 151–161 (2002).
- Siebenhaar, F. et al. The search for mast cell and basophil models - Are we getting closer to pathophysiological relevance? *Allergy Eur. J. Allergy Clin. Immunol.* 70, 1–5 (2015).
- Ackerman, S.J. To be, or not to be, an eosinophil: that is the ??? *Blood* 122, 621–623 (2016).
- Stevens, W. W., Kim, T. S., Pujanauski, L. M., Hao, X. & Braciale, T. J. Detection and quantitation of eosinophils in the murine respiratory tract by flow cytometry. *J. Immunol. Methods* 327, 63–74 (2007).
- Stöger, J. L., Goossens, P. & de Winther, M. P. J. Macrophage heterogeneity: relevance and functional implications in atherosclerosis. *Curr. Vasc. Pharmacol.* 8, 233–48 (2010).
- Stritt, S. et al. Rap1-GTP-interacting adaptor molecule (RIAM) is dispensable for platelet integrin activation and function in mice. *Blood* 125, 219–222 (2015).
- Teodosio, C. et al. The immunophenotype of mast cells and its utility in the diagnostic work-up of systemic mastocytosis. *J. Leukoc. Biol.* 97, 49–59 (2015).
- Torrero, M. N., Larson, D., Hübner, M. P. & Mitre, E. CD200R surface expression as a marker of murine basophil activation. *Clin. Exp. Allergy* 39, 361–369 (2009).
- Travers, J. & Rothenberg, M. E. Eosinophils in mucosal immune responses. *Mucosal Immunol.* 8, 464–75 (2015).
- Tung, J. W. et al. *Modern Flow Cytometry: A Practical Approach.* Clin. Lab. Med. 27, 453–468 (2007).
- Tung, J. W., Parks, D. R., Moore, W. A., Herzenberg, L. A. & Herzenberg, L. A. New approaches to fluorescence compensation and visualization of FACS data. *Clin. Immunol.* 110, 277–283 (2004).
- Turner, J. D. et al. Circulating CD14^{bright}CD16⁺ 'Intermediate' Monocytes Exhibit Enhanced Parasite Pattern Recognition in Human Helminth Infection. *PLoS Negl. Trop. Dis.* 8, e2817 (2014).

- Vermaelen, K. & Pauwels, R. Accurate and simple discrimination of mouse pulmonary dendritic cell and macrophage populations by flow cytometry: Methodology and new insights. *Cytom. Part A* 61, 170–177 (2004).
- Platelets and Megakaryocytes. (Springer New York, 2012). doi: 10.1007/978-1-61779-307-3
- Stem Cells and Aging. (Springer New York, 2013). doi: 10.1007/978-1-62703-317-6
- Eosinophils. (Springer New York, 2014). doi: 10.1007/978-1-4939-1016-8
- Wang, X. S., Yip, K. H., Sam, S. W. & Lau, H. Y. A. Buffy coat preparation is a convenient source of progenitors for culturing mature human mast cells. *J. Immunol. Methods* 309, 69–74 (2006).
- Wang, X., Sze Wing Sam, Kwok Ho Yip & Lau, H. Y. A. Functional characterization of human mast cells cultured from adult peripheral blood. *Int. Immunopharmacol.* 6, 839–847 (2006).
- Wernersson, S. & Pejler, G. Mast cell secretory granules: armed for battle. *Nat. Rev. Immunol.* 14, 478–94 (2014).
- Wilcox, D. A. Megakaryocyte and megakaryocyte precursor related gene therapies. *Blood* 127, 1260–1269 (2016).
- Winter, O. et al. Megakaryocytes constitute a functional component of a plasma cell niche in the bone marrow. *Blood* 116, 1867–1875 (2010).
- Wong, K. L. et al. The three human monocyte subsets: Implications for health and disease. *Immunol. Res.* 53, 41–57 (2012).
- Wright, G. J. et al. Characterization of the CD200 Receptor Family in Mice and Humans and Their Interactions with CD200. *J. Immunol.* 171, 3034–3046 (2003).
- Wynn, T. a, Chawla, A. & Pollard, J. W. Macrophage biology in development, homeostasis and disease. *Nature* 496, 445–55 (2013).
- Xue, J. et al. Transcriptome-Based Network Analysis Reveals a Spectrum Model of Human Macrophage Activation. *Immunity* 40, 274–288 (2014).
- Yallapu, M. M., Jaggi, M. & Chauhan, S. C. Curcumin nanoformulations: a future nanomedicine for cancer. *Drug Discov. Today* 17, 71–80 (2012).
- Yang, Q. & Bhandoola, A. The development of adult innate lymphoid cells. *Curr. Opin. Immunol.* 39, 114–120 (2016).
- Yeaman, M. R. Platelets: at the nexus of antimicrobial defence. *Nat. Rev. Microbiol.* 12, 426–37 (2014).
- Zaynagetdinov, R. et al. Identification of myeloid cell subsets in murine lungs using flow cytometry. *Am. J. Respir. Cell Mol. Biol.* 49, 180–189 (2013).

Zhang, J., Socolovsky, M., Gross, A. W. & Lodish, H. F. Role of Ras signaling in erythroid differentiation of mouse fetal liver cells: functional analysis by a flow cytometry – based novel culture system. *Differentiation* 102, 3938–3946 (2003).

Ziegler-Heitbrock, L. Reprint of: Monocyte subsets in man and other species. *Cell. Immunol.* 291, 11–15 (2014).

Zimring, J. C. et al. A novel mouse model of red blood cell storage and posttransfusion in vivo survival. *Transfusion* 49, 1546–1553 (2009).