

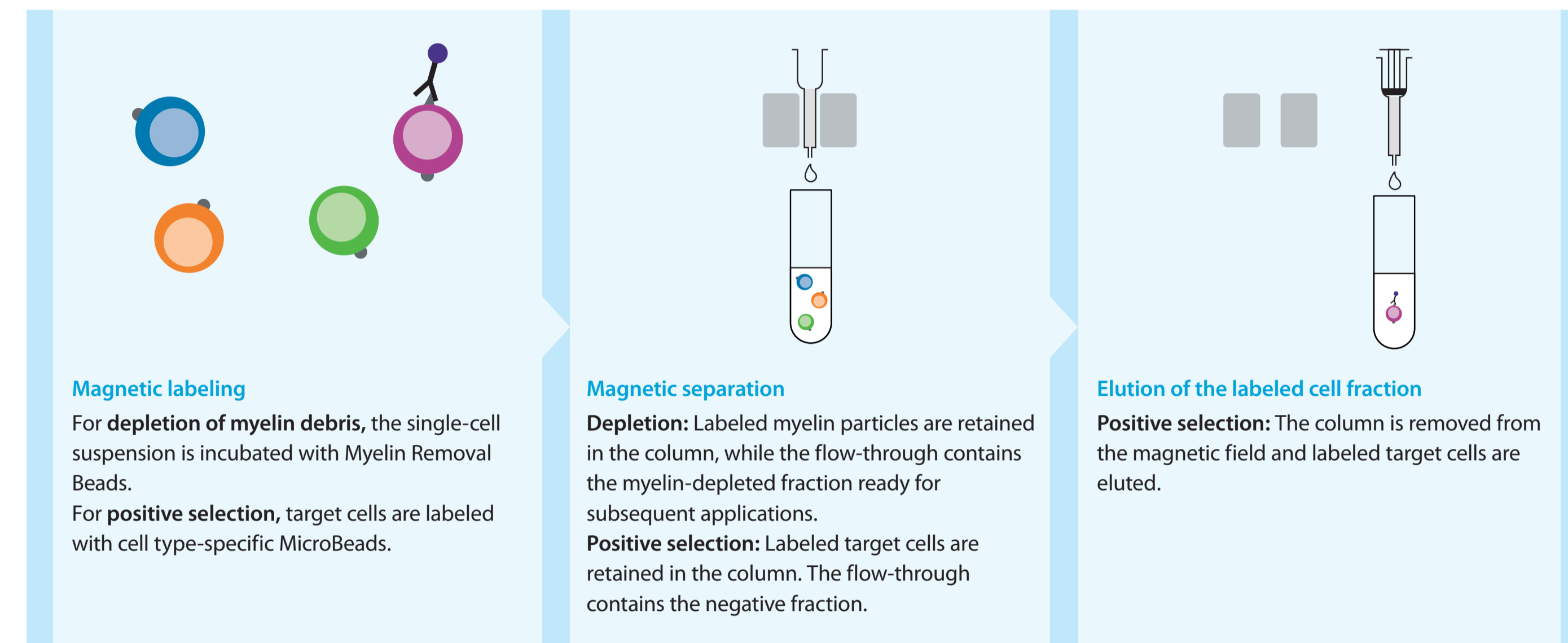
Removal of myelin debris after tissue dissociation optimizes immunomagnetic sorting of cells from postnatal brain

Sandy Reiß, Ina Herzig, Andreas Bosio, Sandra Pennartz | Miltenyi Biotec GmbH, Bergisch Gladbach, Germany

Introduction

Myelin, a specialized membrane, ensheathes and insulates axons in the peripheral and central nervous system. In mice and rats, myelination begins around birth in the spinal cord and is completed in the brain during the first postnatal month¹. In humans, myelin formation starts during the second half of fetal life in the spinal cord, peaks during the first year post-natally and can continue until 20 years of age².

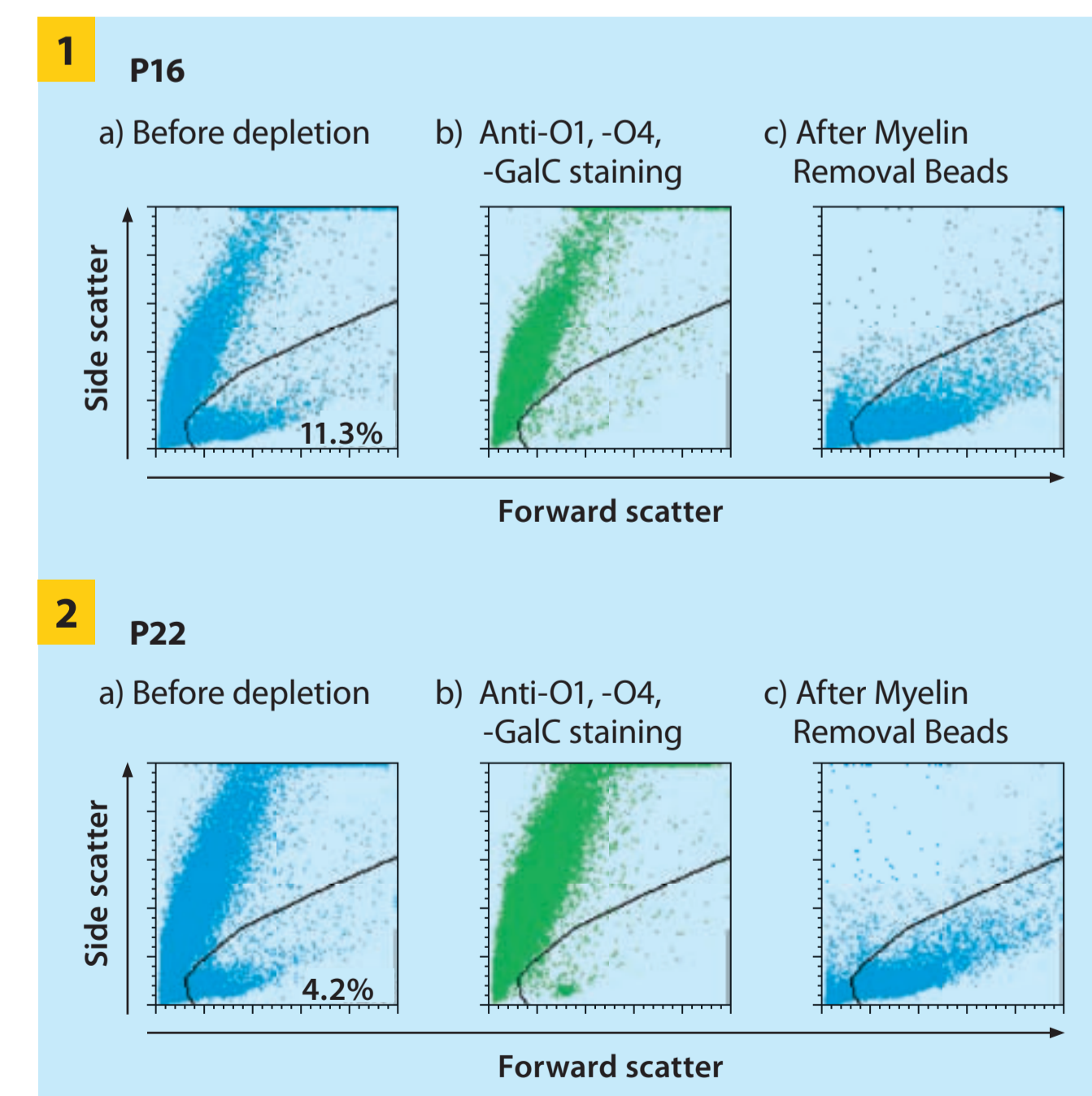
When myelin containing neural tissue is dissociated, large quantities of myelin debris are generated. Myelin debris in single-cell suspensions was reported to considerably impair cell isolation and antibody staining^{3,4}. So far, sucrose^{5,6} has been used for the elimination of myelin. Here, we present a solution for myelin depletion based on magnetic cell sorting, MACS[®] Technology (see illustration below).



MACS[®] Technology for cell separation. Either certain cells—or components—can be depleted, or cells can be positively selected for presence of a certain surface molecule.

Results

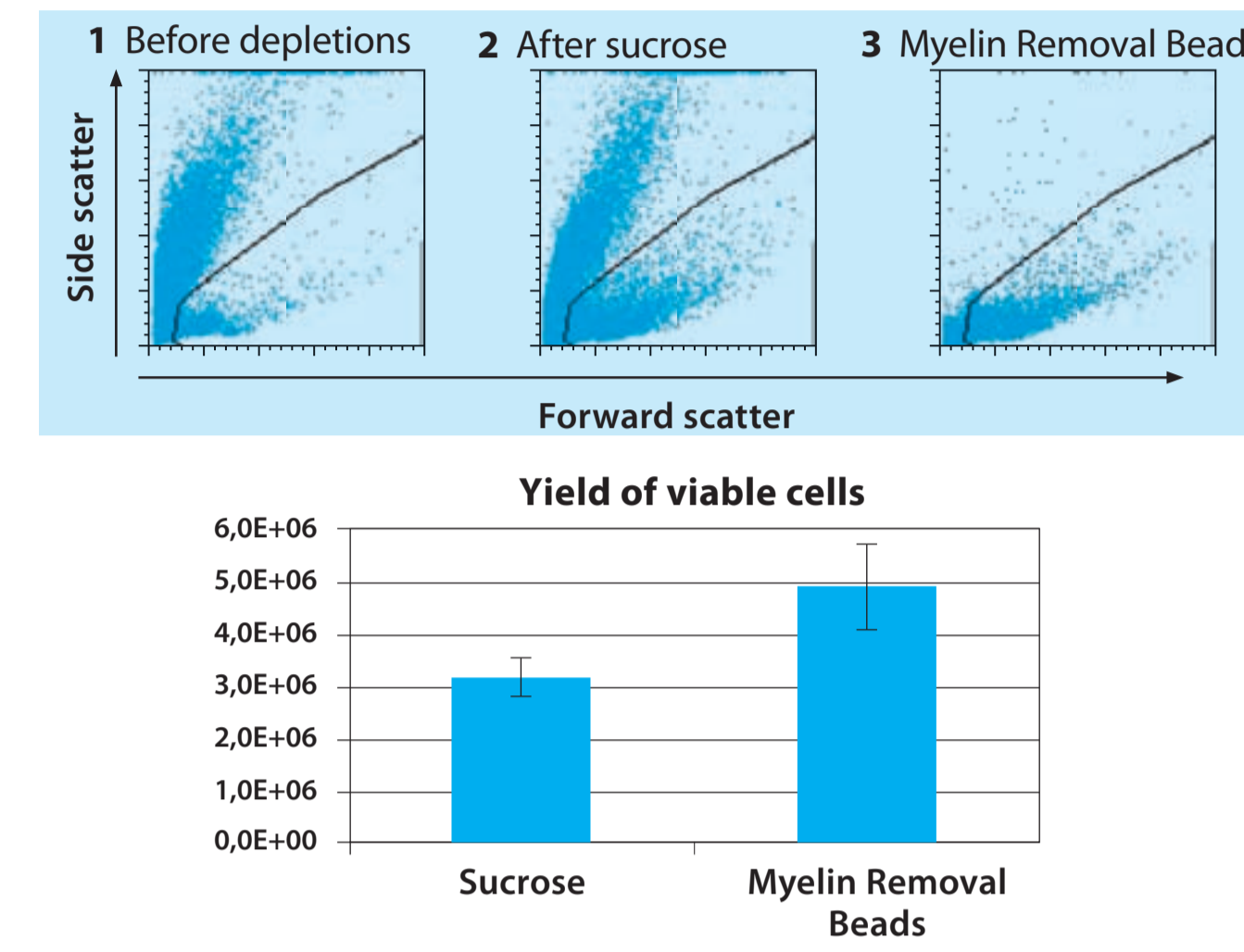
1 Myelin debris is abundant in single-cell suspensions



Mouse brains were dissociated using the Neural Tissue Dissociation Kit-P (NTDK-P, Miltenyi Biotec). Flow cytometrical analysis of the resulting single-cell suspensions shows the distribution of cells and myelin debris (a): forward scatter (cell size) and side scatter (granularity).

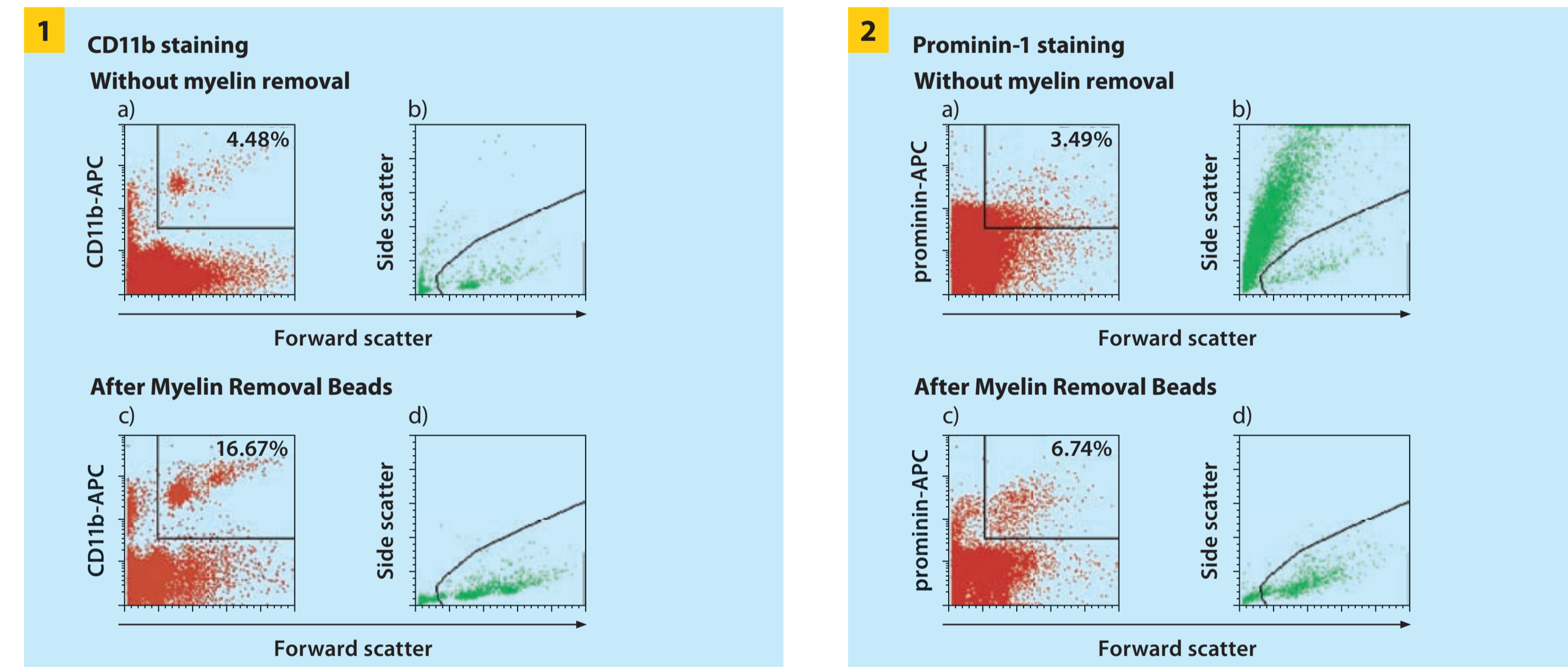
Single-cell suspensions from postnatal day (P) P16 (1) and P22 (2) mouse brains consist of large amounts of myelin membrane fragments, other cellular debris and only 11% and 4% cells, respectively. Particles outside the selected regions are positive for anti-O1, anti-O4 and anti-GalC antibodies that recognize glycolipids specifically contained in myelin membranes (b). Myelin Removal Beads used with LD Columns and VarioMACS[™] Separators efficiently remove myelin debris (c).

2 Myelin Removal Beads are more efficient than sucrose solution for myelin debris removal



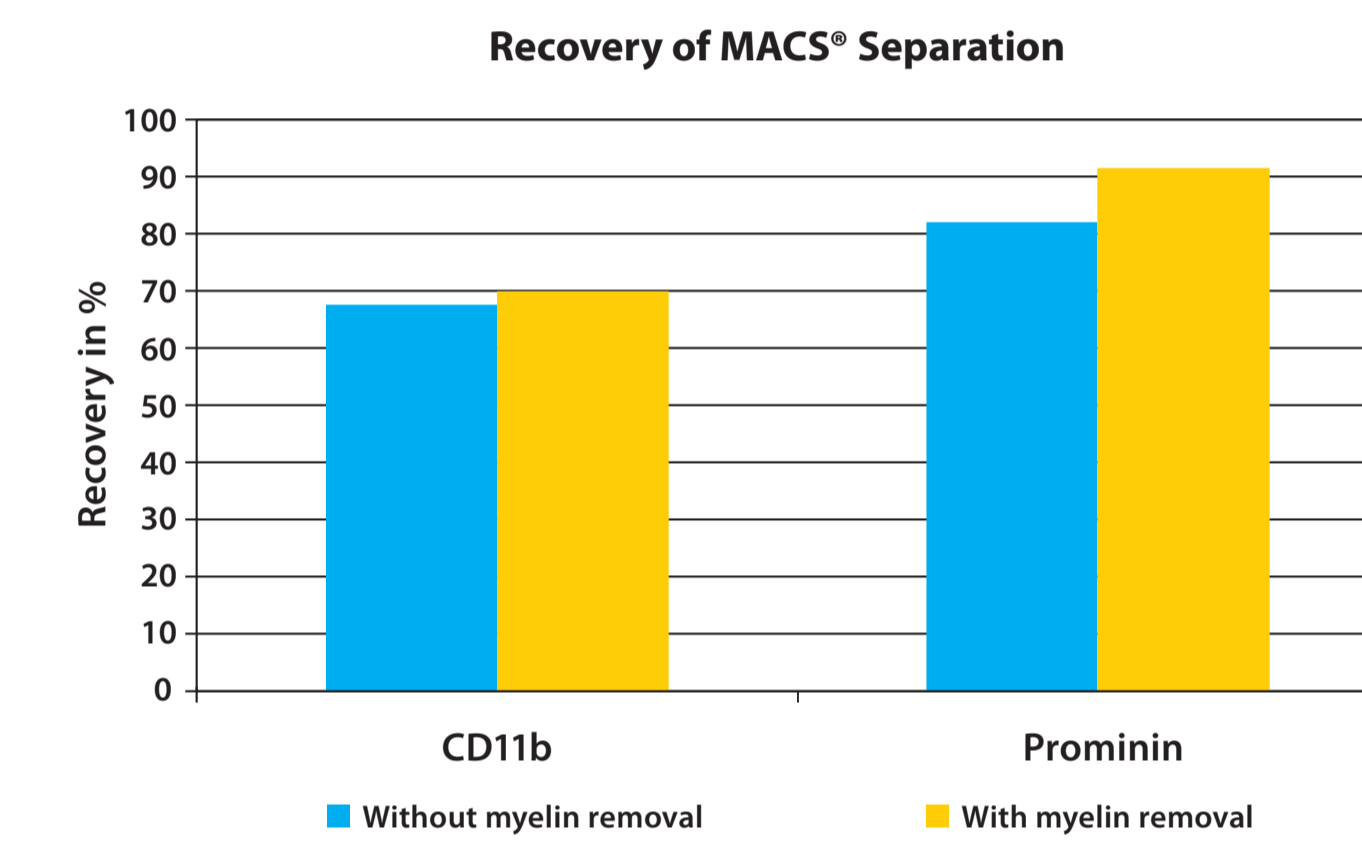
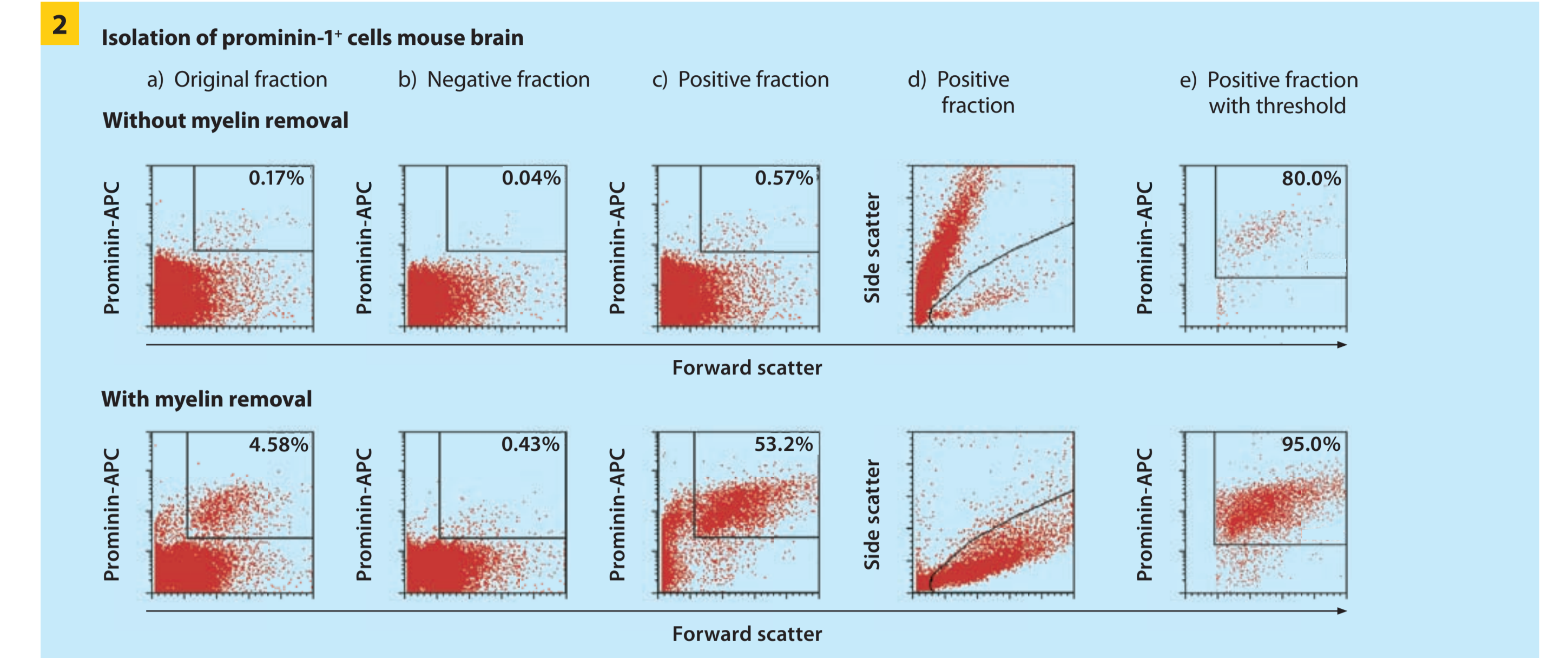
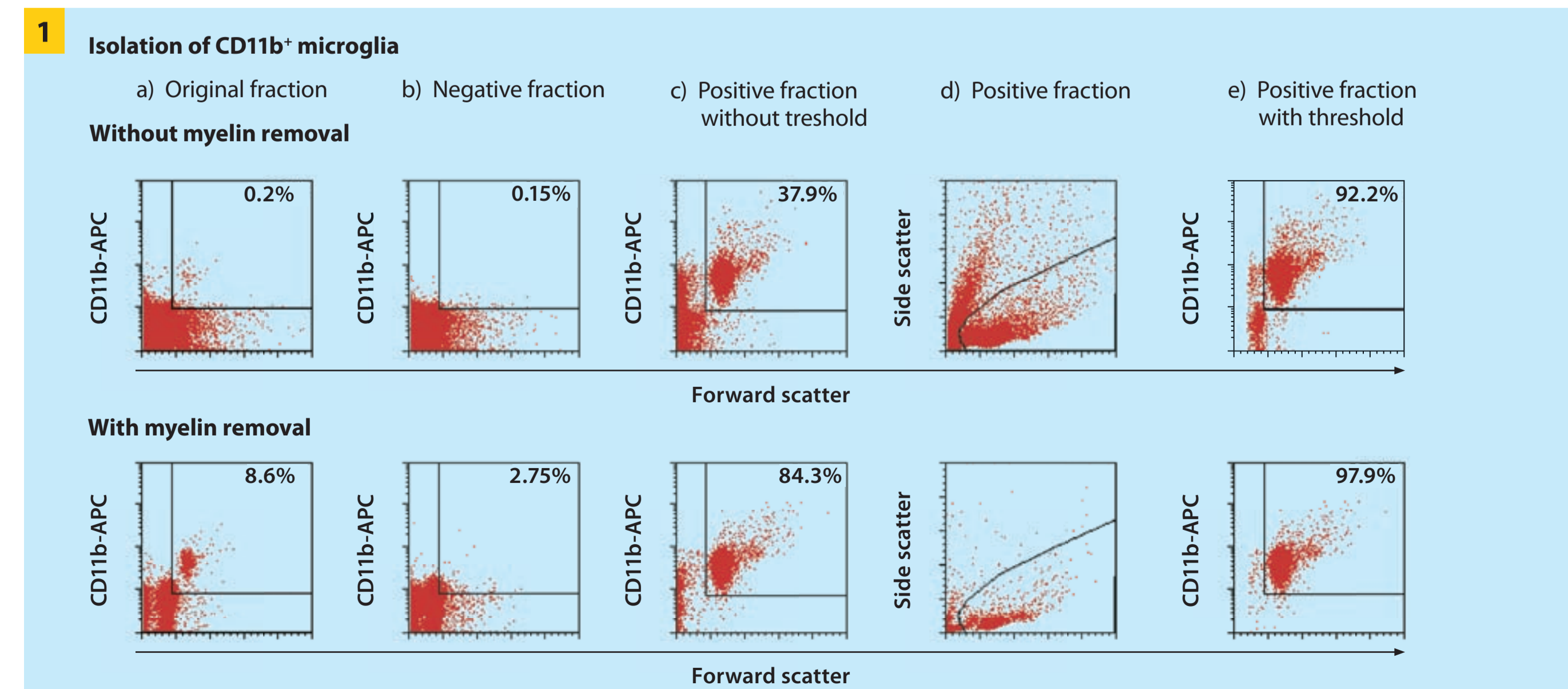
Myelin Removal Beads (Miltenyi Biotec) deplete myelin debris more thoroughly than sucrose solution. Dot plots show the amount of myelin debris and cells in forward and side scatter before depletion (1), after treatment with 10 mL of 0.9 M sucrose and centrifugation at 850×g for 10 minutes⁶ (2) and after depletion using Myelin Removal Beads (3). Dead cells were labeled with PI and are not shown in the dot plots (n = 5). The diagram illustrates that Myelin Removal Beads produce a higher recovery of viable cells from a starting population of 1×10⁷ cells than treatment with sucrose solution.

3 Removal of myelin debris improves antibody stainings



1×10⁶ cells from a single-cell suspension of P22 mouse brain were stained with CD11b-APC (1) or anti-Prominin-1-APC (2) without and with previous myelin removal using Myelin Removal Beads. Dot plots 1a+c and 2a+c show that in samples with previous myelin removal, higher percentages of CD11b- and Prominin-1-positive cells are stained, respectively. Dead cells were excluded using PI. Only the positive cells along with positive debris are displayed in dot plots 1b+d and 2b+d in side and forward scatter.

4 Removal of myelin debris increases efficiency of MACS[®] Separations



For MACS Separations using CD11b-MicroBeads (1a–e) or Anti-Prominin-1 MicroBeads (2a–e), P18 or P22 mouse brain, respectively, was dissociated using the NTDK-T. 1×10⁷ cells from a single-cell suspension were directly used for separation or were submitted to myelin depletion using Myelin Removal Beads. Comparing the separations from samples without and with myelin removal demonstrates that the recovery is higher for samples with previous myelin removal. These data demonstrate a higher efficiency of cell separations in the absence of myelin debris.

Conclusions

- Myelin Removal Beads efficiently deplete myelin debris
- The recovery of viable cells after Myelin Removal Beads is higher than after sucrose solution
- Removal of myelin significantly improves the results of immunostainings and cell isolation experiments

References

- 1 Simons, M. and Trotter, J. (2007) Wrapping it up: the cell biology of myelination. *Curr. Opin. Neurobiol.* 17: 533–540.
- 2 Baumann, N. and Pham-Dinh, D. (2001) Biology of oligodendrocyte and myelin in the mammalian central nervous system. *Physiol. Rev.* 81: 871–927.
- 3 Pfenninger, C. V., Roschupkina, T., Hertwig, F., Kottwitz, D., Englund, E., Bengzon, J., Jacobsen, S. E., and Nuber, U. A. (2007) CD133 is not present on neurogenic astrocytes in the adult subventricular zone, but on embryonic neural stem cells, ependymal cells, and glioblastoma cells. *Cancer Res.* 67: 5727–5736.
- 4 Tham, C. S., Lin, F. F., Rao, T. S., Yu, N., and Webb, M. (2003) Microglial activation state and lysophospholipid acid receptor expression. *Int. J. Dev. Neurosci.* 21: 431–443.
- 5 Johansson, C. B., Momma, S., Clarke, D. L., Risling, M., Lendahl, U., and Frisen, J. (1999) Identification of a neural stem cell in the adult mammalian central nervous system. *Cell* 96: 25–34.
- 6 Suzchet, S., Arnason, B. G., and Polak, P. E. (1980) Separation of ovine oligodendrocytes into two distinct bands on a linear sucrose gradient. *J. Neurosci. Methods* 3: 7–19.