

A Novel Ultraviolet Dye for Cell Proliferation Analysis in Flow Cytometry

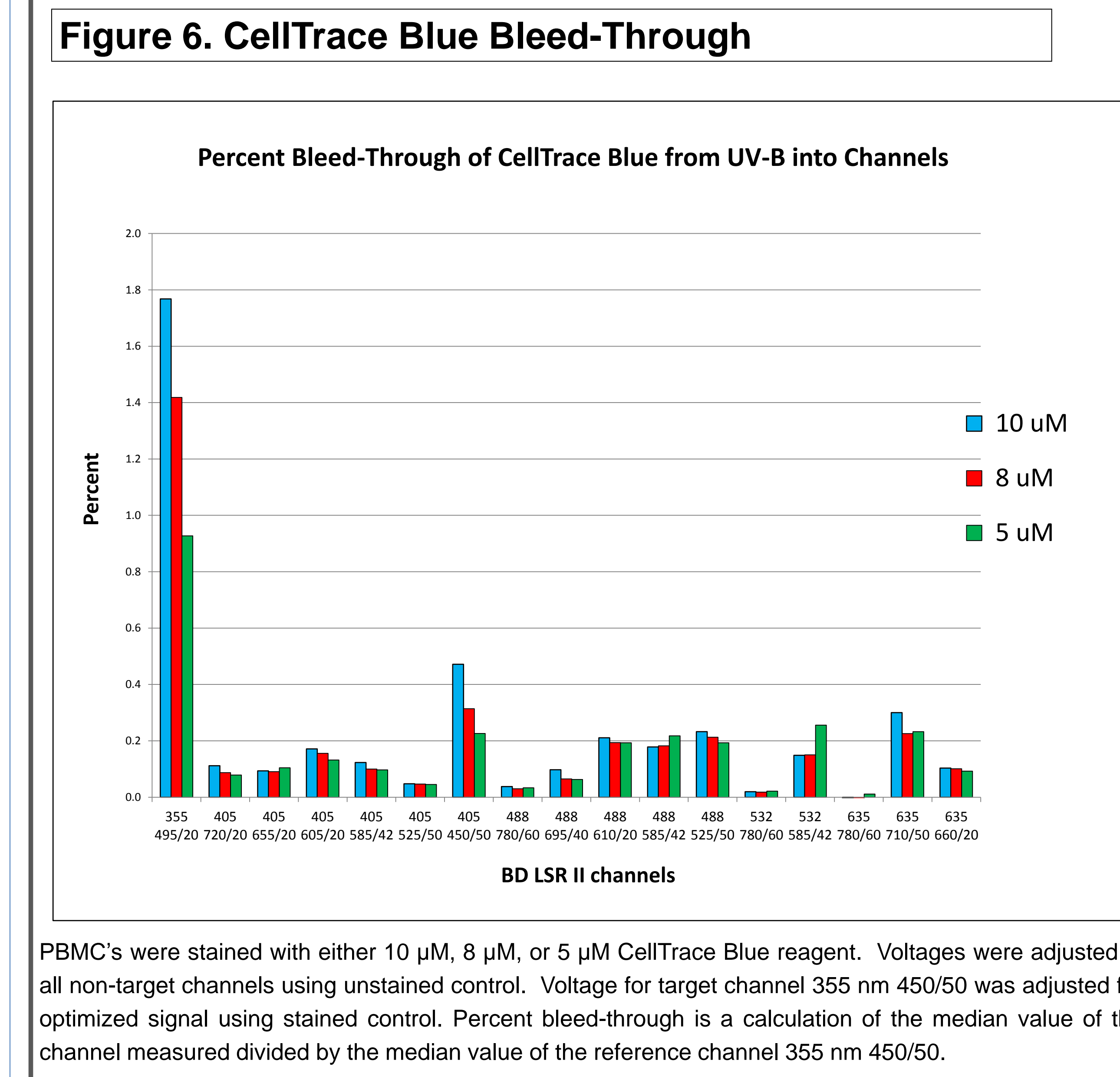
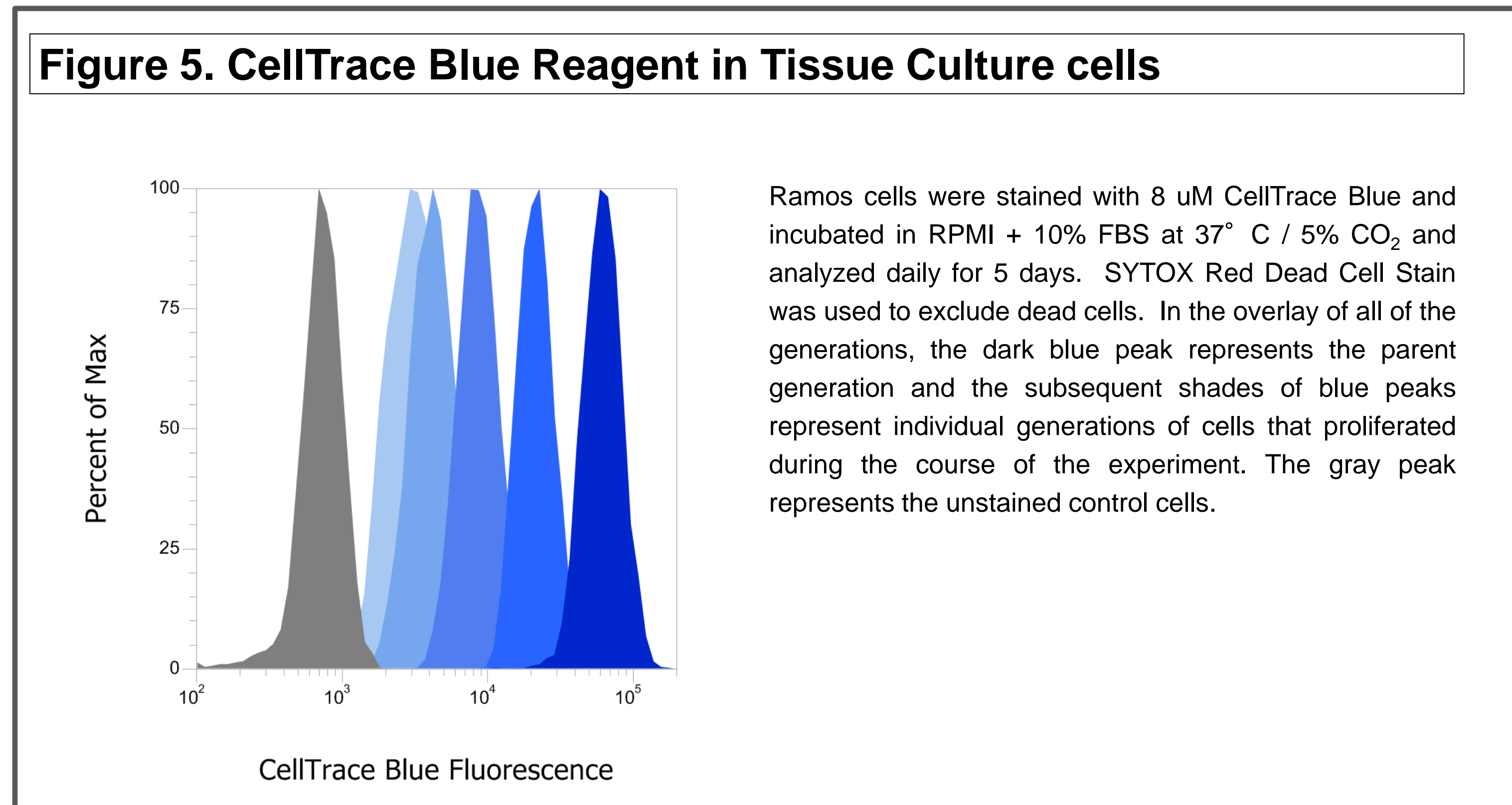
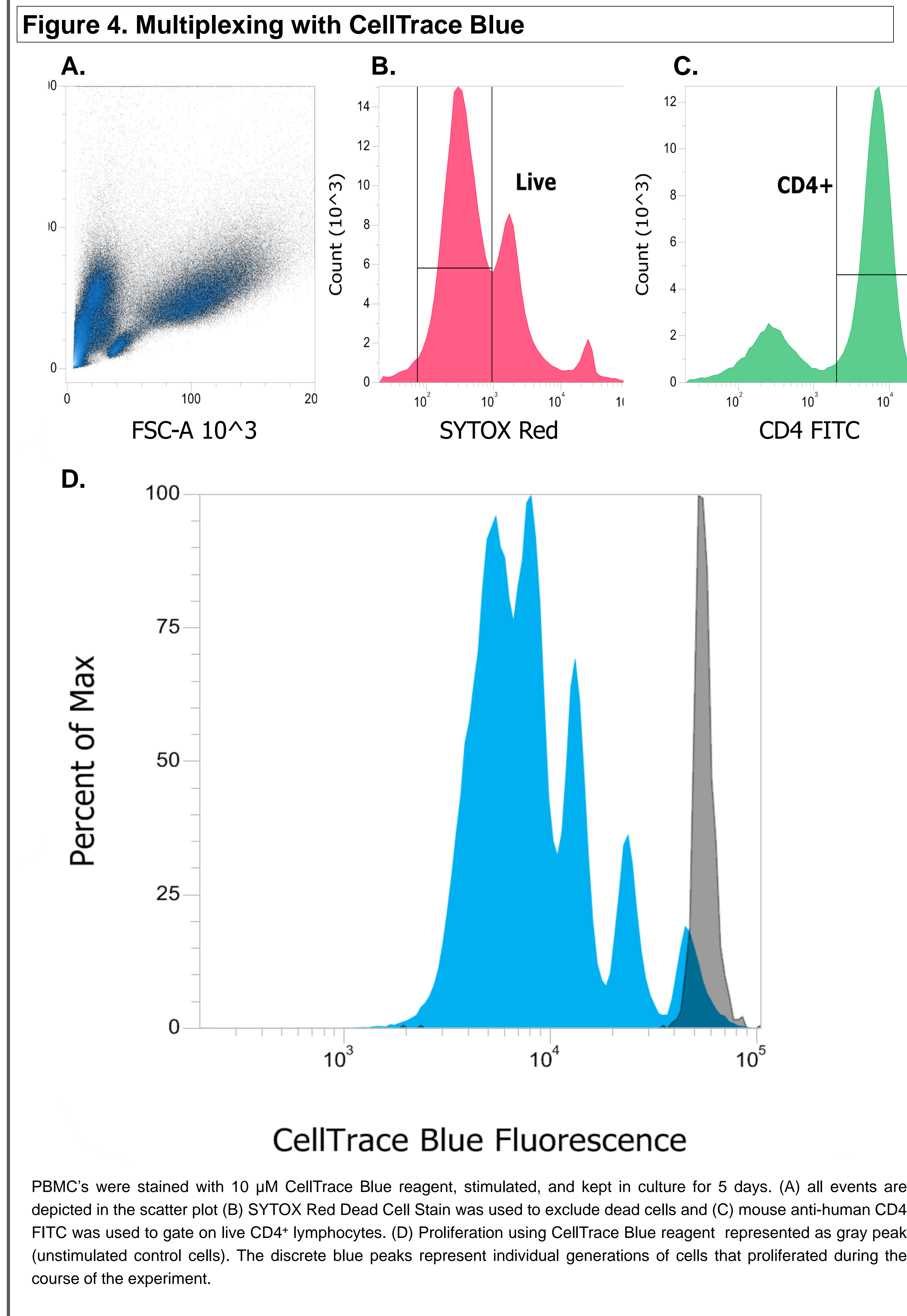
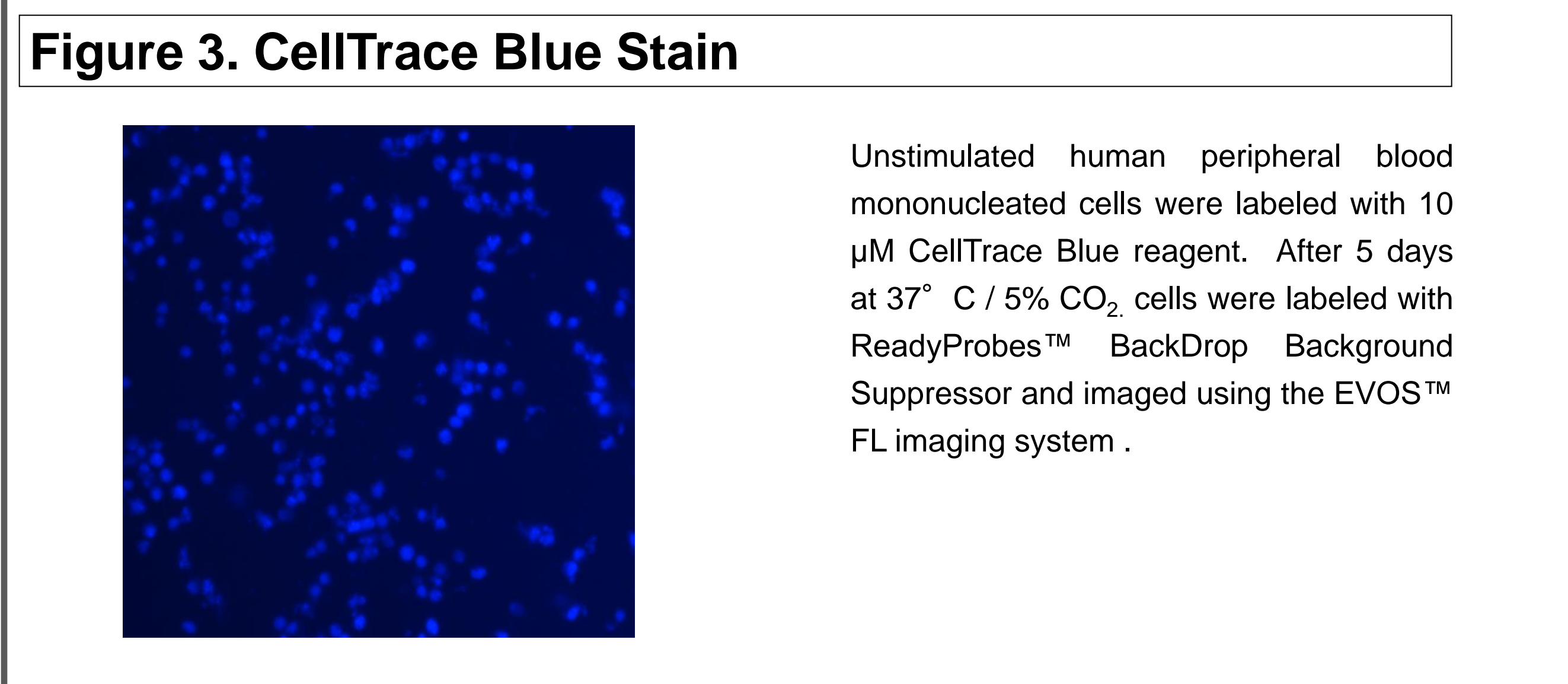
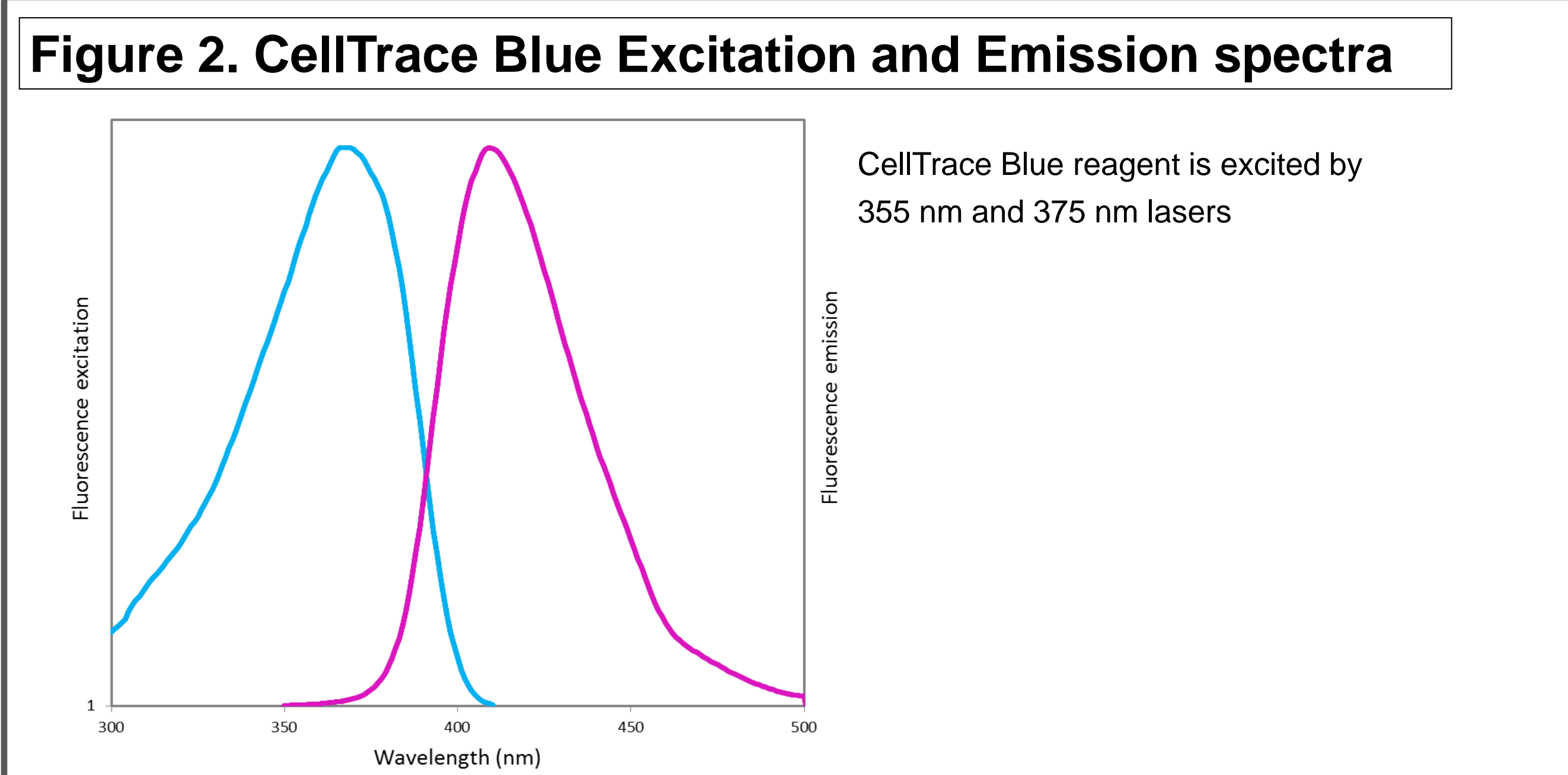
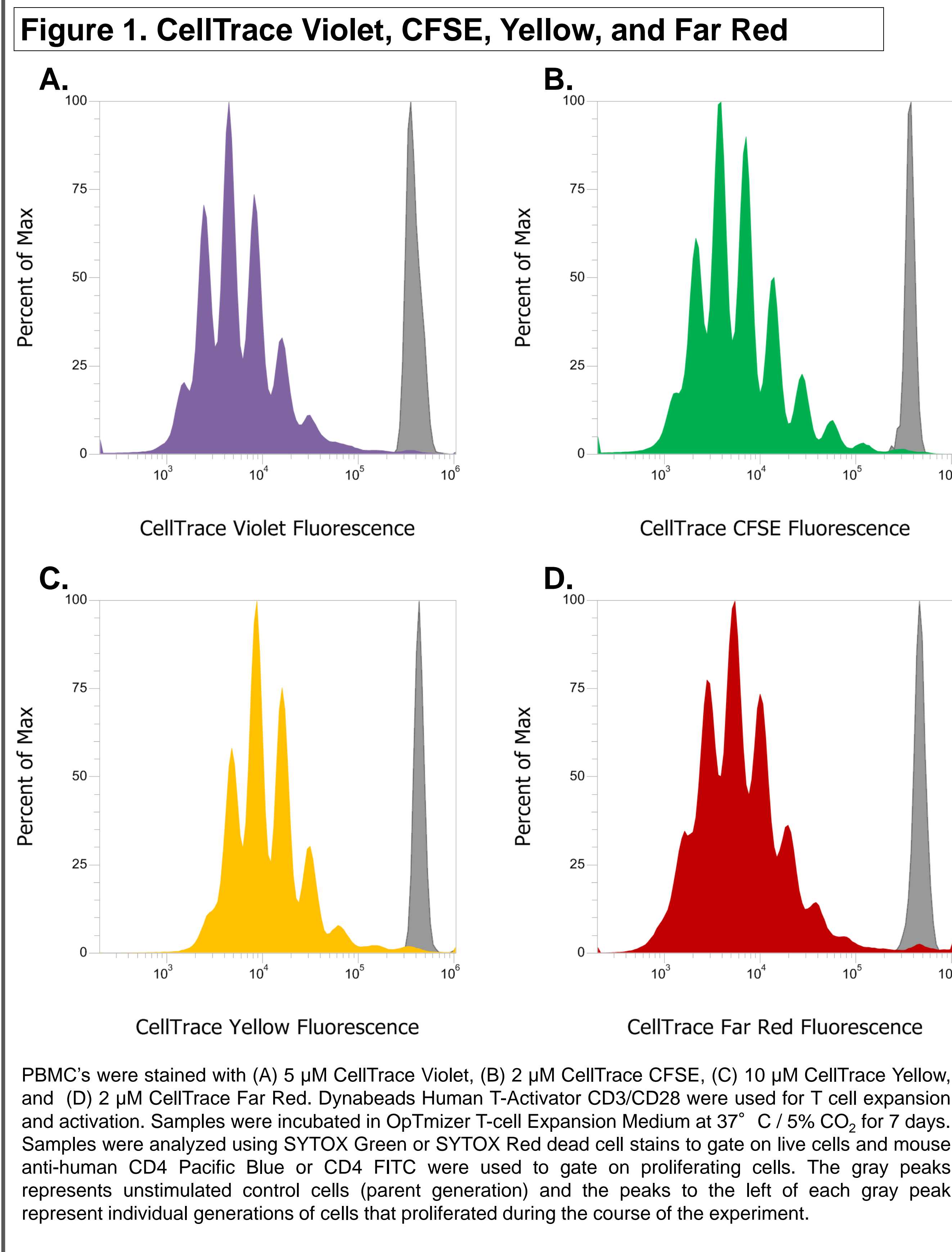
Penny Melquist, Kyle Gee, and Aimei Chen
Thermo Fisher Scientific, 29851 Willow Creek Road, Eugene, OR, USA, 97402

INTRODUCTION

Cell proliferation reagents are crucial for determining cell growth and differentiation and are often used to evaluate compound toxicity and/or inhibition of tumor cell growth. Tools for measuring cell proliferation include probes for analyzing the average DNA content and cellular metabolism in a population, as well as single-cell indicators of DNA synthesis and cell cycle-specific proteins, and tracking population doubling by dye dilution. Successful proliferation analysis by dye dilution requires an extremely bright dye to distinguish fluorescently labeled cells after several cell divisions, however many of the existing reagents require compensation. The CellTrace™ Blue ultraviolet dye frees up channels for other reagents that can be designed into a multi-color panel. The fluorescent staining provided by CellTrace Blue enables the visualization of five generations of proliferating cells. CellTrace Blue covalently binds intracellular proteins where the stable, well-retained fluorescent dye offers a consistent signal, even after several days in cell culture. Division of a cell labeled with a CellTrace Blue results in partitioning of dye between daughter cells and approximately half of the fluorescence intensity in the progeny. When analyzed by flow cytometry using either a 355 nm or 375 nm laser, this partitioning of the CellTrace Blue dye provides a direct indication of cell proliferation without compromising cell health. In this experiment, we will look at how CellTrace Blue complements the existing CellTrace portfolio and demonstrate that the spectral characteristics permit multiplexing with many common flow reagents without compensation, distinct peak determination upon cell proliferation, and performance with tissue culture cells.

MATERIALS AND METHODS

Human peripheral blood mononuclear cells were isolated from whole blood using a Ficoll-Paque Plus density gradient. Cells were stained with 10 μ M CellTrace Blue reagent and resuspended in CTS™ OpTmizer™ T-Cell Expansion media. Stained cells were stimulated to proliferate with 50 μ L CD3/CD28 Dynabeads™ per milliliter cells and incubated at 37° C and 5% CO₂ for 5 days. Cells were removed from culture, washed, and labeled with 5 μ L mouse anti-human CD4 FITC for 20 minutes. Finally, cells were washed and labeled with SYTOX™ Red Dead Cell Stain and data was acquired on the BD LSR II using 355 nm 450/50. Software for the Invitrogen Attune™ NxT was used to analyze the data.



CONCLUSIONS

Results indicate that CellTrace Blue reagent can be used to successfully track multiple generations of proliferating cells, including tissue culture cells. The bright, staining produces minimal fluorescence variation between cells in a population, resulting in the ability to visualize resolve distinct generations by flow cytometry. The ultraviolet excitation at 355 nm or 375 nm make it ideal for multiplexing due to the limited spectral overlap with other common dyes. Results of this experiment suggest that CellTrace Blue reagent is a useful tool to track and characterize cells as they proliferate through several generations.

REFERENCES

- Wallace, et al. *Cytometry Part A* 73A: 1019-1034, 2008.
- Lyons AB, Parish CR. *Journal of Immunological Methods* 1994; 171:131-37.
- Givan AL, et al. *Journal of Immunological Methods* 1999; 230:99-112.

SPECIAL THANKS

Special thanks to Mike O'Grady, Helen Fleisig, Suzanne Buck, Jolene Bradford, Veronica Calderon, Jenny Rosenberger, Hiram Rasmussen, Terrie Goodman, Marcy Wickett, Yi Zhen Hu, & Alece Barley.

TRADEMARKS/LICENSING

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.
Ficoll-Paque is a trademark of GE Healthcare
For research use only. Not for use in diagnostic procedures.

