

An Optimized Fluorescent 9-plex Spatial Proteomics Approach with Spectral Unmixing to Identify Cell Types in the Brain

Denise Niell, Nancie Mooney, Kristi Hamilton, Oggie Golub
Thermo Fisher Scientific, Eugene, Oregon, United States

Abstract

The ability to uniquely identify cell types increases greatly with the number of biomarkers that can be used concurrently. Traditional immunohistochemistry (IHC) approaches have limitations in terms of resolving spectrally distinct signals for multiple targets. We have developed a system for flexible labeling and amplification of protein targets for fluorescent IHC with spectral unmixing, which enables the separation of overlapping spectral signatures and the accurate identification of eight biomarkers within a single tissue section. Using the EVOS S1000 Spatial Imaging System to generate high-resolution images of whole tissue sections combined with quantitative image analysis software, we can measure the spatial distribution of distinct cell populations based on identified biomarkers. This allows for the specific localization of neuronal and glial subpopulations within a single tissue section. These results highlight the advantages of using a spectrally unmixed 9-plex sample analysis (8 biomarkers and a nuclear counterstain) to investigate the spatial relationships between cell types in the brain.

Multiplex Labeling

Aluora™ Spatial Amplification Kits

Aluora kits use tyramide-like signal amplification labeling technique for enhanced sensitivity and intensity. Combining the brightness of Invitrogen™ fluorescent dyes with poly-HRP-mediated signal amplification labeling permits high-fidelity multiplexing of a variety of validated antibody clones.

Materials and methods

Sample Preparation

Formalin-fixed, paraffin-embedded (FFPE) mouse coronal brain tissue sections (5µm) were obtained from Zyagen Inc. These slides were then processed using a Leica Bond RXm and stained with primary antibodies (table right) and the Aluora Spatial Amplification kits, as described in Figure 1. Images were acquired and spectrally unmixed on the EVOS S1000.

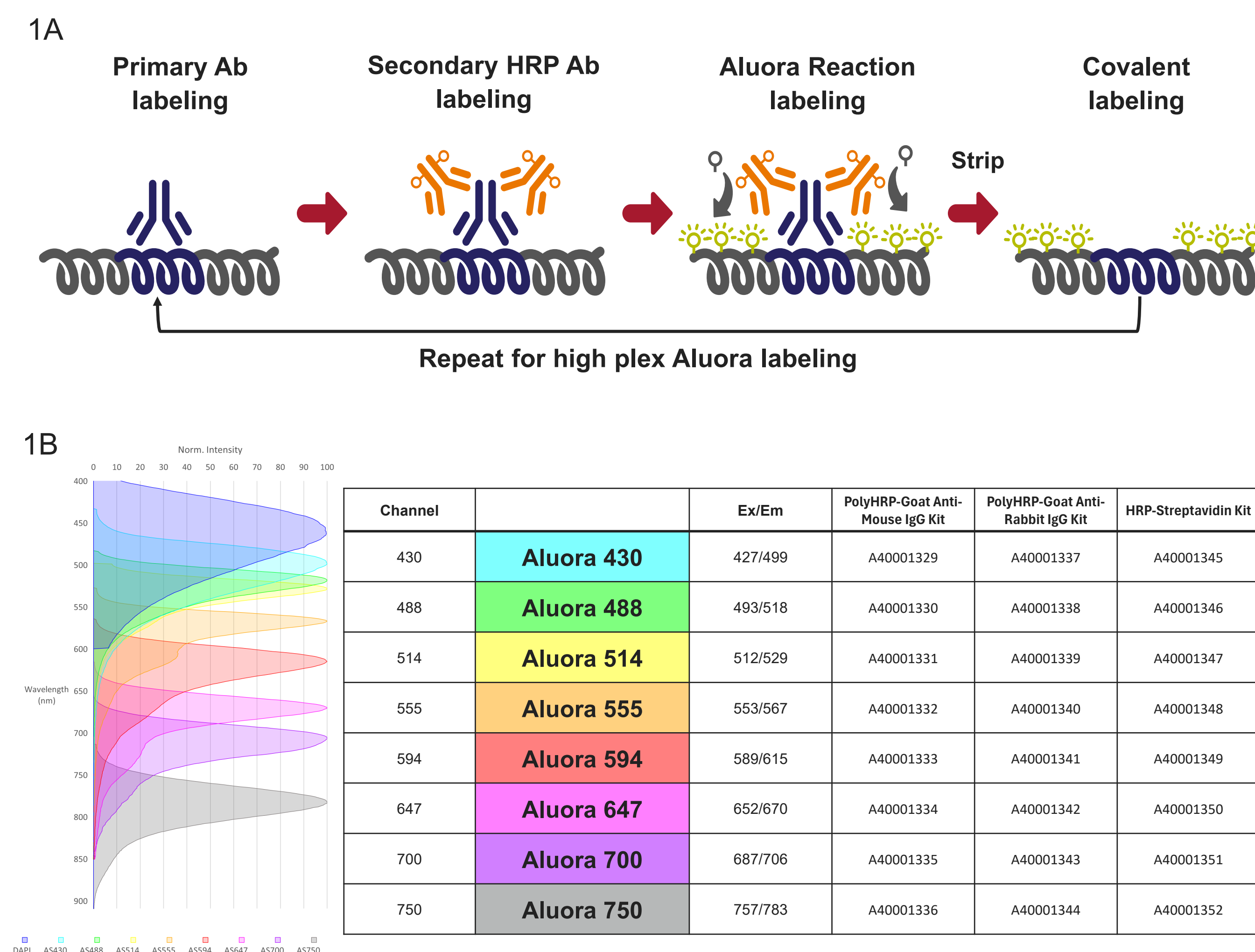
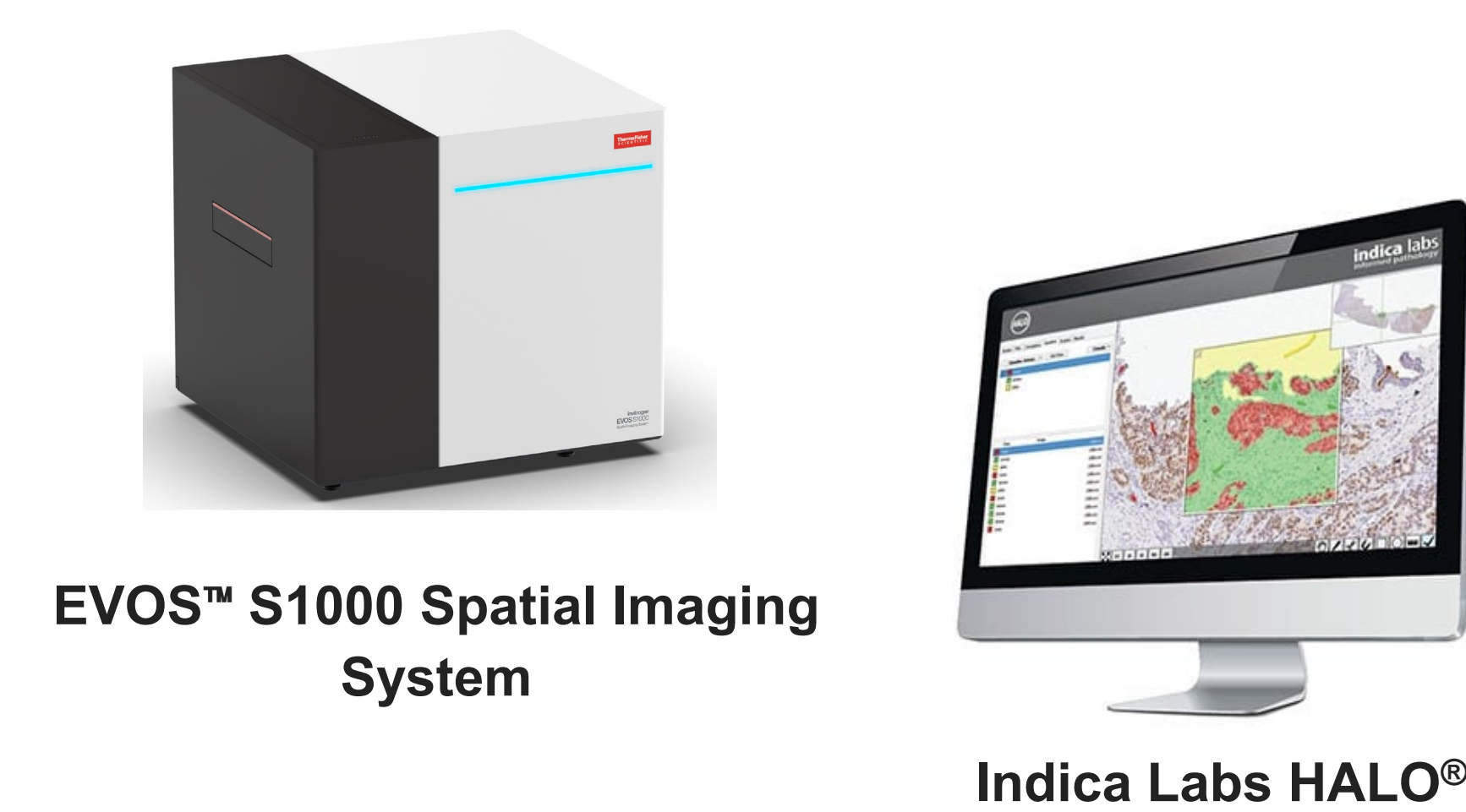


Figure 1. Multiplex labeling workflow with the Aluora kits uses horseradish peroxidase (HRP) for enzymatic conversion of the Aluora dyes to covalently bind tissue on and surrounding the protein epitope targeted primary antibodies. A) Diagram of workflow. B) Emission spectra and table of Aluora products.



Data Analysis

Analysis of the multiplex immunofluorescence stitched image was performed on the Indica Labs HALO® (version 4.0.5107.318) software. 81,926 cells were identified across a coronal section using the Indica Labs-HighPlex FL version 4.2.14, and the Halo AI Nuclei Seg V2-FL classifier.

Learn more at thermofisher.com/spatial-biology

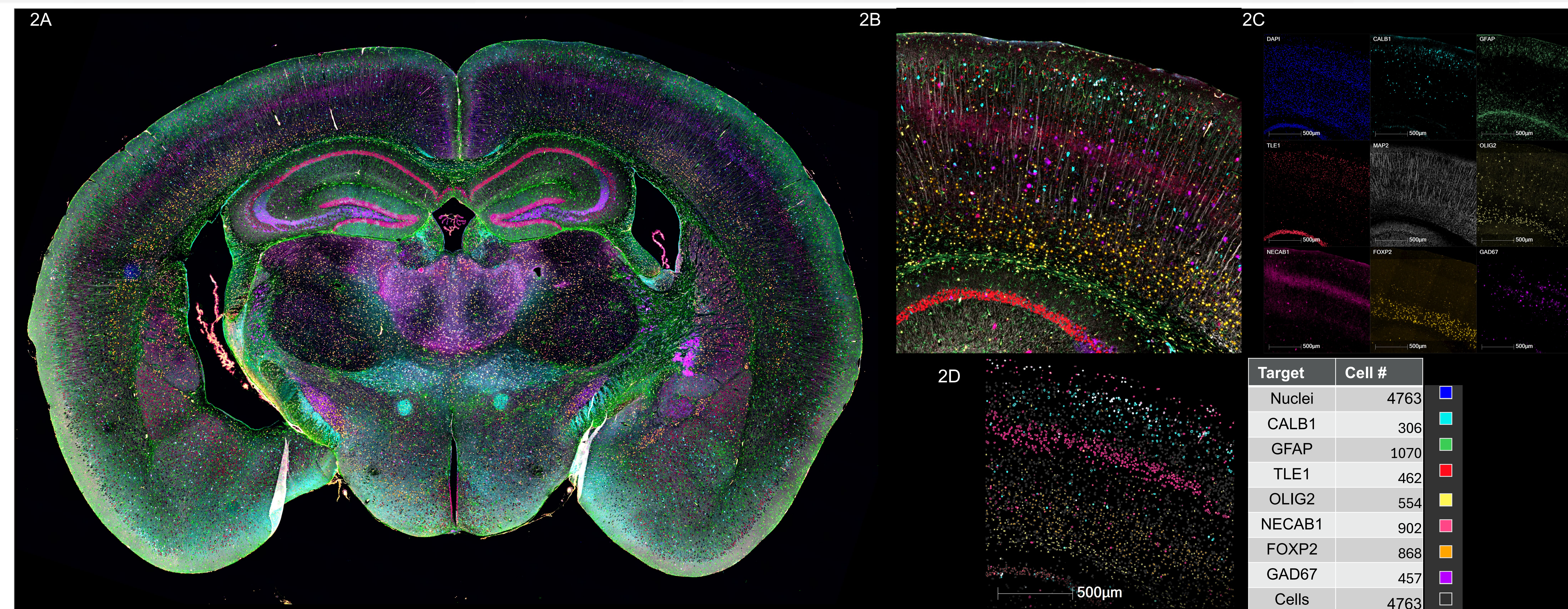


Figure 2. Mouse brain tissue stained with the 8-plex Aluora spatial amplification assay and DAPI. Mouse coronal brain tissue section was processed and stained with the Aluora spatial amplification 8-plex assay, 20x Imaging and spectral unmixing performed on the EVOS S1000 Spatial Imaging System. A) Composite image, B) zoomed-in inset, C) individual unmixed channels, D) Quantification of target positive cells identified using HALO®.

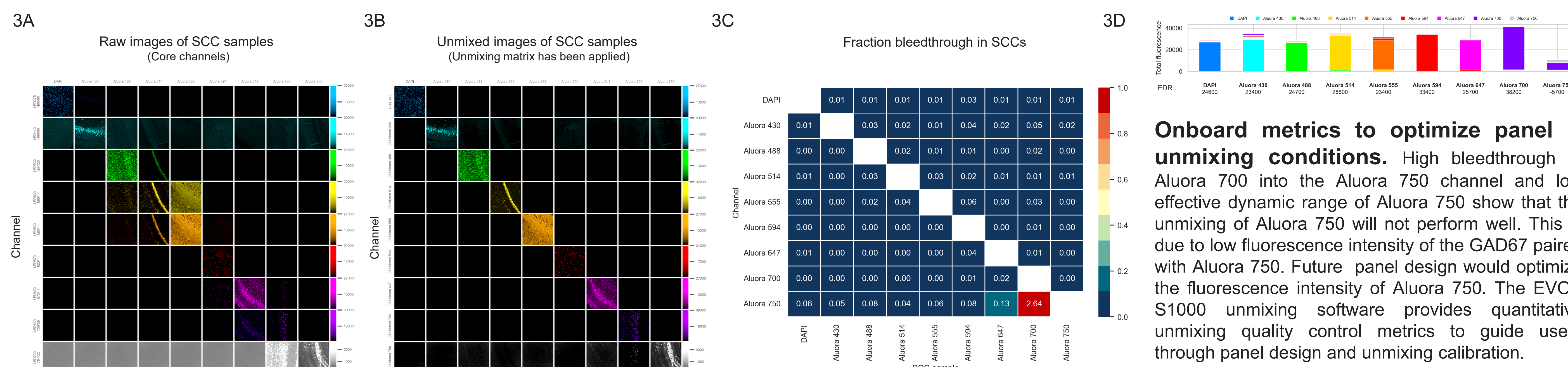


Figure 3. Spectral unmixing enables high-plex spatial biology imaging. A) Single color control (SCC) samples showing spectral overlap in raw images where fluorophore is detected by neighboring core channels. B) Application of the unmixing matrix separates signals into their respective channels. C) Quantification of fluorescence in “Unmixed images”. Fraction bleed-through is the ratio of the signal observed in off target channels to the signal of the target fluorophore after unmixing has occurred. D) Effective dynamic range (EDR) = Target fluorescence – bleedthrough.

Conclusions

We developed a highly flexible and sensitive Aluora labeling method capable of 8-rounds of protein labeling. Combined with the EVOS S1000, which enables spectral unmixing of overlapping dyes, we demonstrate the ability to distinguish the unique staining patterns of 8 different biomarkers and DAPI. Onboard metrics assist with the optimization of complex panel design. These high-resolution images allow for precise quantification of 8 targets and identification of cell types in an entire mouse brain section at 20x magnification.

Trademarks/licensing

© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. For Research Use Only. Not for use in diagnostic procedures.