

Antibody Development and Validation in the Era of Spatial Proteomics

biotechne® /
R&D SYSTEMS™

Ashley Oliver, Kristine Trueman, Jodi Hagen, Alex Kalyuzhny, and Nidhi Nath
R&D Systems, a Bio-Techne brand, Antibody Development, Minneapolis, MN

Learn more about
multiplex IF (mIF)
antibody validation.



01 // Overview

- Spatial proteomics can create high resolution map of large number of proteins at subcellular resolution within tissues, enabling understanding of complex cellular processes.
- Visualizing large array of proteins within complex tissue environment requires highly sensitive and specific antibodies.
- In collaboration with Lunaphore, we have deployed processes to meet the exacting standards for antibodies on the spatial biology COMET™ platform.
- Antibodies designed for purpose will allow broad adoption of spatial proteomics and multiomics in research, drug discovery and diagnostic fields.

02 // COMET™ an Automated Sequential Immuno-fluorescence (seqIF™) Platform for Spatial Biology

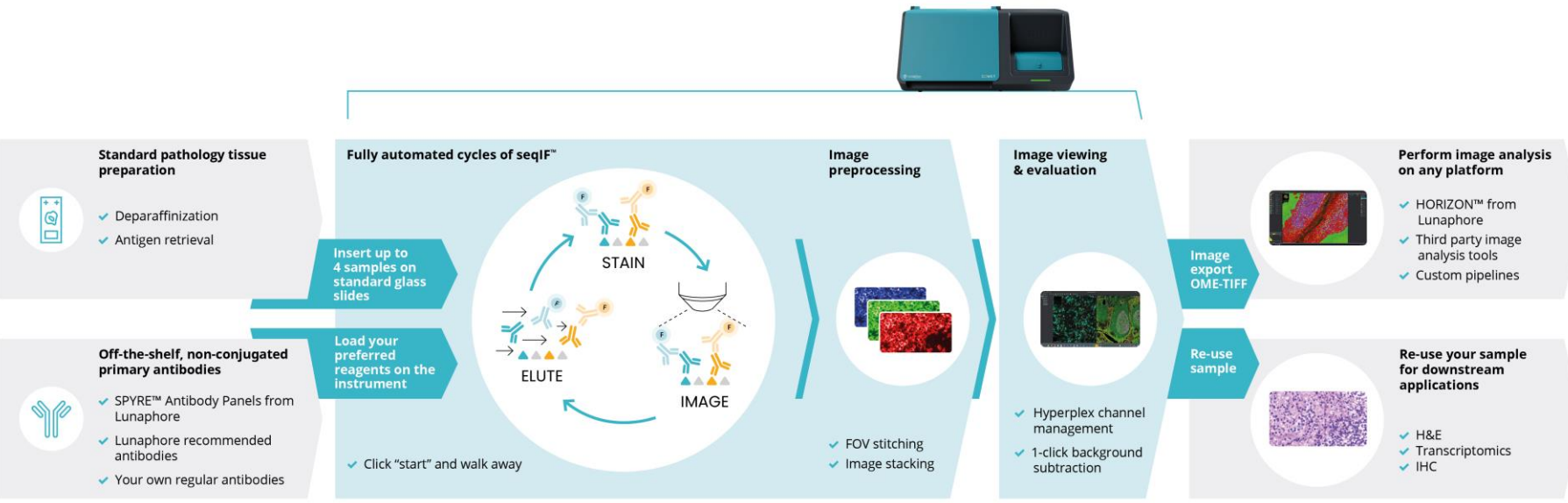


Figure 1. COMET™ - automated staining & imaging platform for multiplex spatial biology

- Unmatched multiplex throughput with walk-away automation
- Rapid and flexible panel development using label-free antibodies
- True reproducibility and tissue preservation

03 // Expanded COMET™ Antibody Discovery Workflow

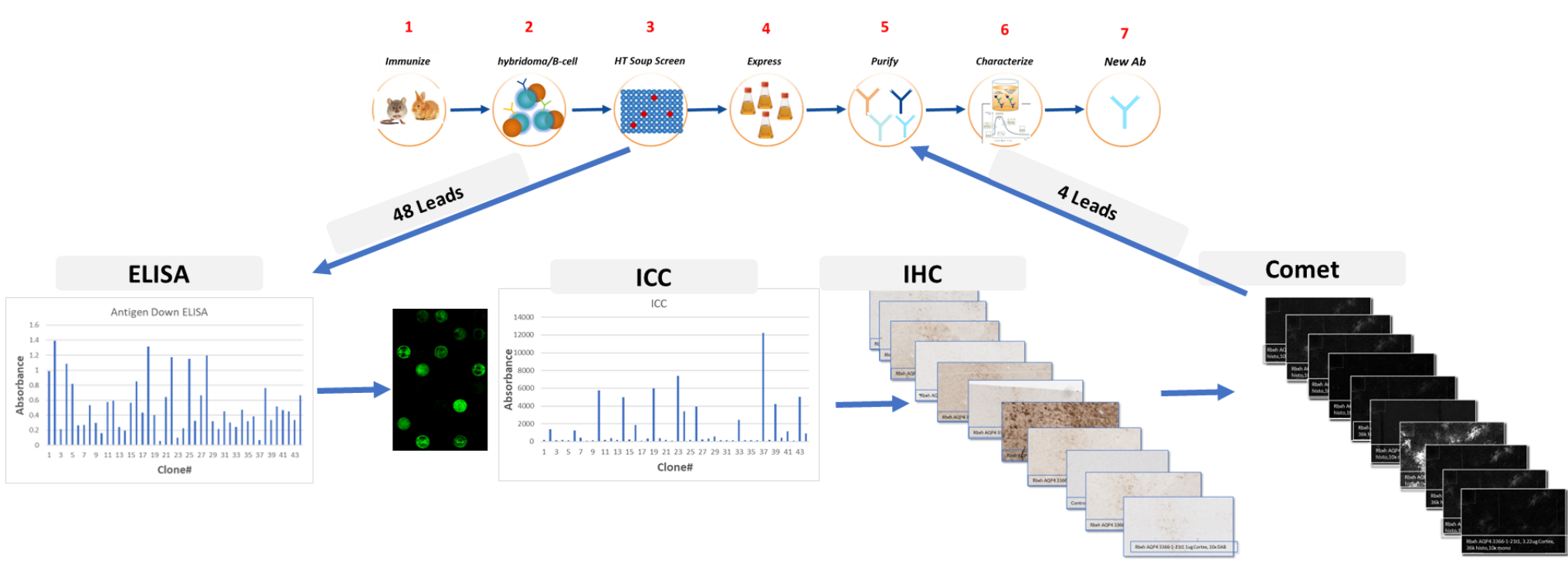


Figure 2. Antibody discovery and validation workflow to select sensitive and specific antibodies for COMET™

- Improved methodology for immunization
- Expanded primary screening to identify lead candidates
- Enhanced testing for platform specific performance

04 // COMET™ Specific Optimization

For best performance on COMET™, antibodies are optimized for, concentration, elution condition, elution cycles, sensitivity, and specificity.

Smooth Muscle Actin (α -SMA) Antibody Optimization

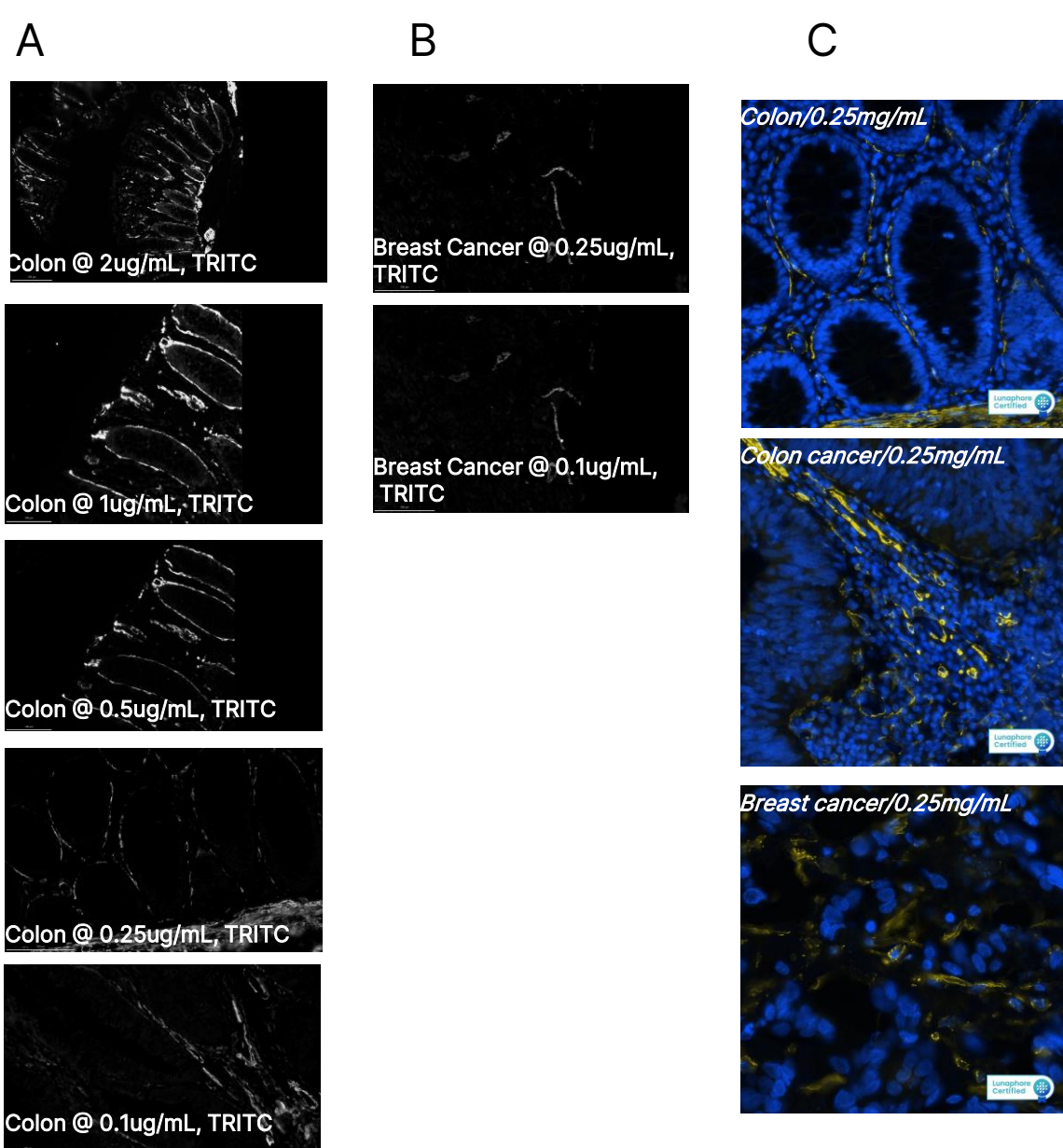


Figure 3. Representative images showing impact of SMA antibody concentration on tissue staining in colon (A) and breast cancer (B). (C) Final images using optimized antibody concentration of 0.25 ug/mL in colon, colon cancer, and breast cancer tissues

05 // COMET™ Antibodies Optimized for Specificity

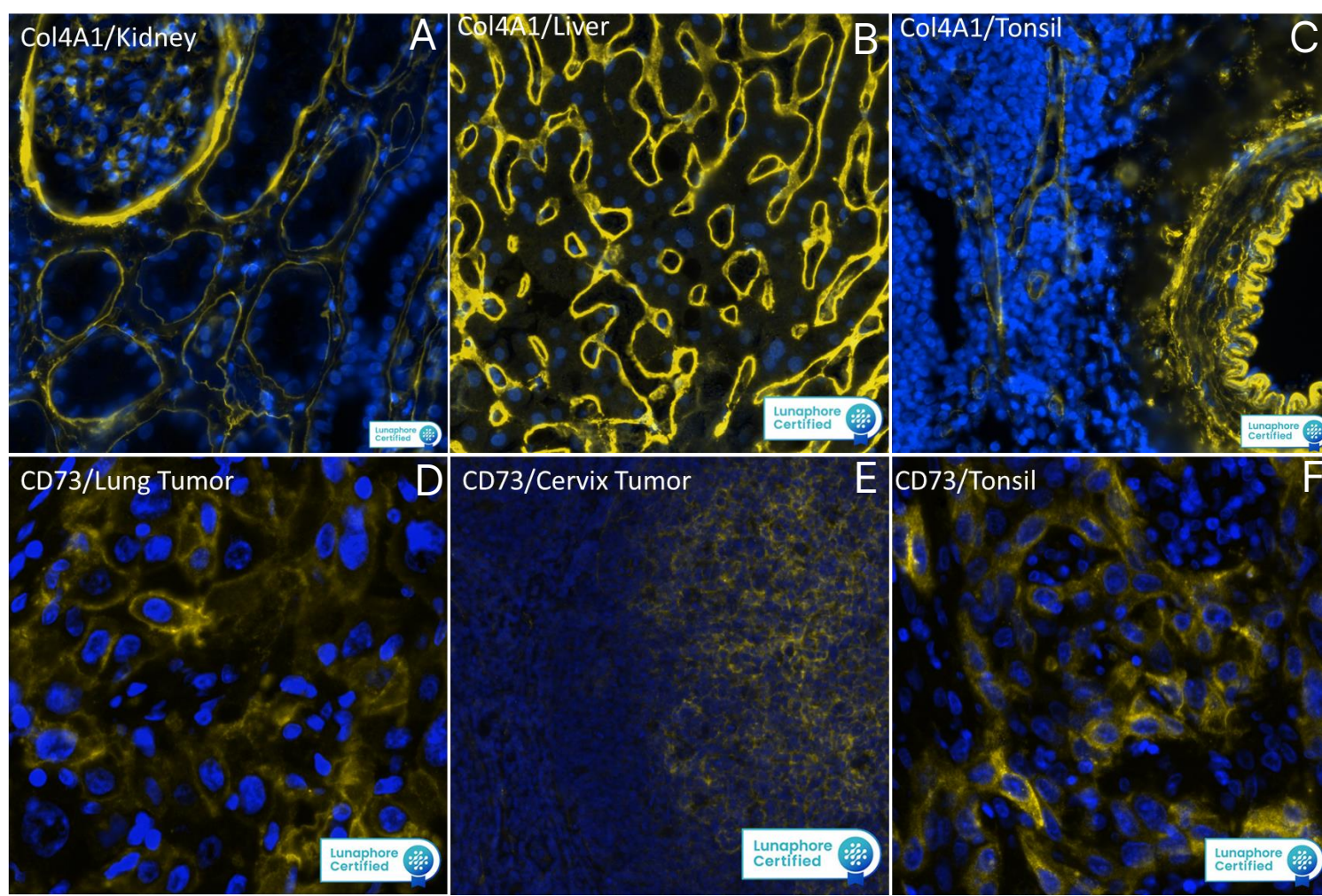


Figure 4. COMET™ optimized antibodies are tested for specificity using multiple tissues. Collagen type IV alpha 1 chain (COL4A1t) localization in basement membrane in kidney (A); liver (B) and tonsil (C). CD73 localization in lung tumor (D); cervix tumor (E); and tonsil (F).

06 // Multiplexed Tissue Image using COMET™ Optimized Antibodies

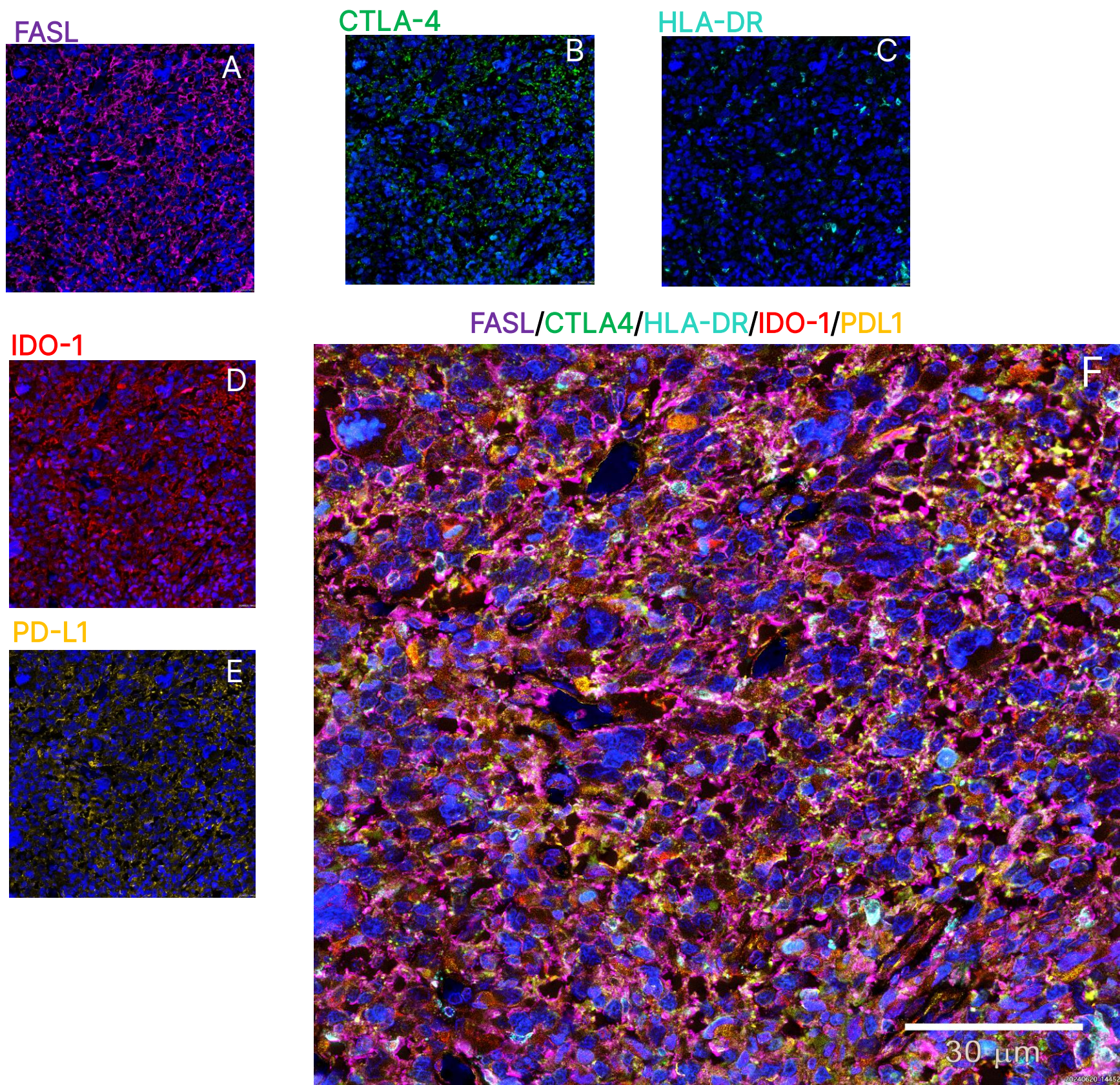


Figure 5. Multiplexed fluorescence imaging in glioblastoma tissues for five different protein markers-FASL (A), CTLA-4 (B), HLA-DR (C), IDO-1 (D) and PD-L1 (E). (F) Composite image highlighting the spatial distribution of five markers on the tissue.

Antibodies used were optimized for COMET™

07 // Results and Conclusions

- We demonstrate that enhanced development and testing strategies can meet the needs for high quality spatial biology antibodies.
- COMET™ platform stands out for multiplexed imaging by eliminating the need for antibody conjugation while enabling automation and high-throughput analysis.
- We are actively expanding and qualifying our antibody catalog to provide customers with an increasingly diverse and robust portfolio of COMET™ tested antibodies for their research needs.

biotechne® // Global Developer, Manufacturer, and Supplier of High-Quality Reagents, Analytical Instruments, and Precision Diagnostics.

INCLUDES R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen® Lunaphore™