



SignalStar™ multiplex immunohistochemistry is a flexible spatial technology with fully validated protocols.

Background

Exploring spatial biology through multiplex immunohistochemistry (mIHC) empowers researchers to delve into the intricate functions, arrangements, and interactions of cells that shape the tumor microenvironment (TME) during disease progression or in response to therapy. SignalStar introduces an innovative mIHC assay employing oligo-conjugated antibodies and a matrix of fluorescent oligonucleotides, amplifying up to 8 targets within a single FFPE tissue.

Methods

Rigorous validation ensures optimal conjugation efficiency and ensures antibody sensitivity and specificity. The myeloid compartment of the TME was assessed using antibodies targeting CD11c, SIRPα, CD163, CD206, CD68, CD45, HLA-DRA, and Pan-Keratin. The simultaneous application of all antibodies is followed by amplifying the signal of four targets in the initial imaging round. Fluorescent signal was removed, and the signal of 4 additional antibodies were amplified in the second round of imaging. SignalStar staining was quantitatively compared to the established chromogenic gold standard to confirm reproducibility across modalities and between replicates.

Results

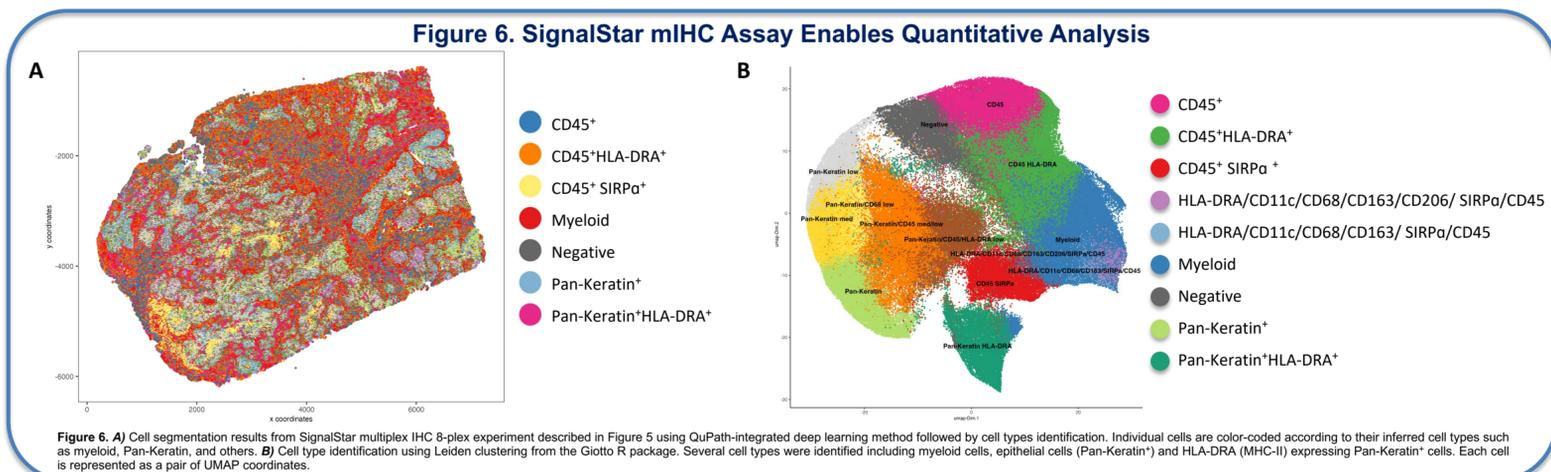
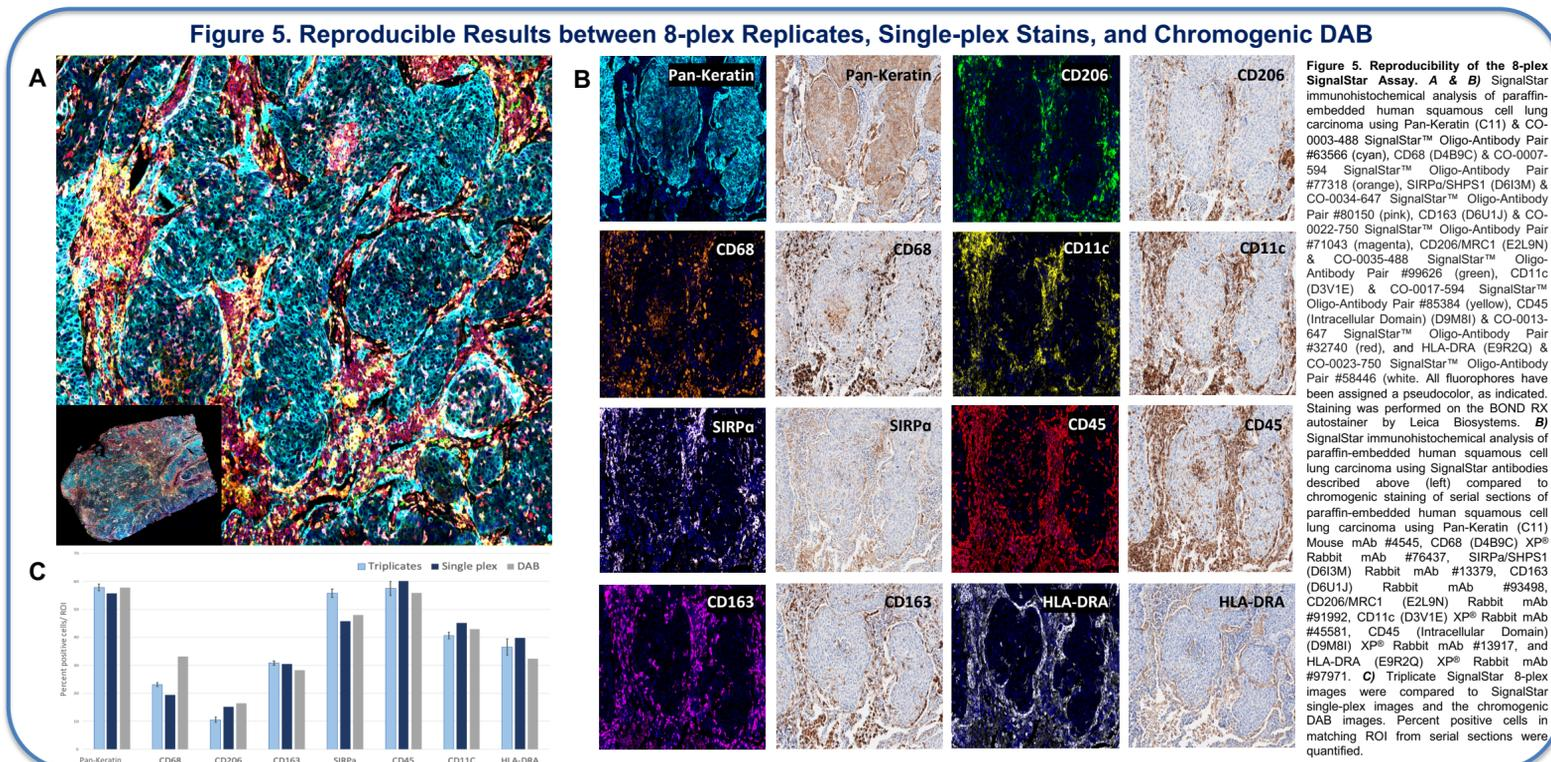
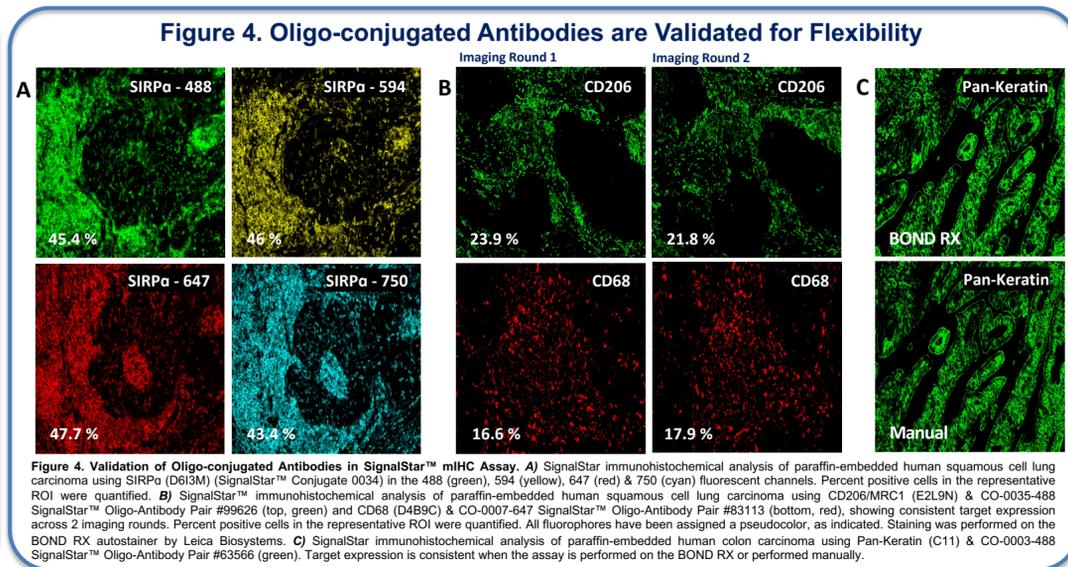
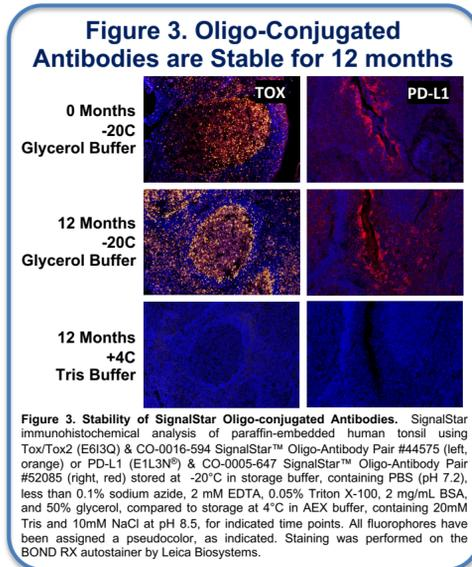
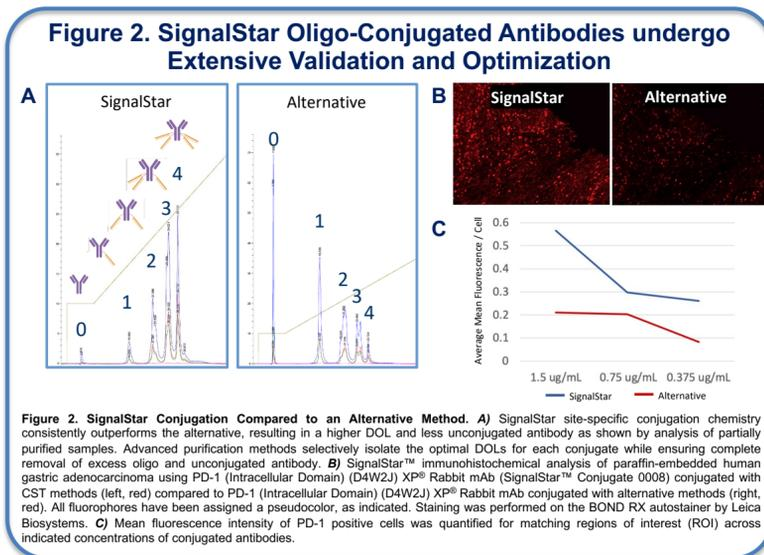
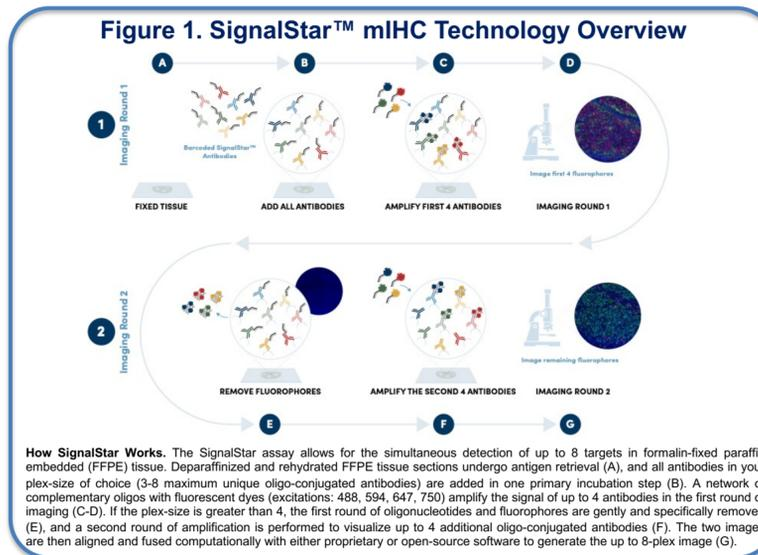
Our findings demonstrate the consistent performance of each antibody across all available fluorescent channels in multiple rounds of imaging regardless of manual or automated protocols, as exemplified by SIRPα and Pan-Keratin. Image alignment and quantification of target signal frequency and co-localization was performed using QuPath. Anticipated cell subsets expressing multiple biomarkers, such as quadruple CD206+CD68+CD163+SIRPα+ cells, were successfully identified. The SignalStar mIHC technology facilitates adaptable panel design without lengthy optimization while upholding precision and reproducibility in quantifiable phenotypic data for comprehensive immune cell interrogation within the TME.

Conclusions

Spatial imaging is increasingly important to understand the different biological aspects of the TME, including the cells present and the organization and function of their biomarkers. This can be critical to understanding the differences between diseased and healthy tissue. Current technology is laborious and time consuming, particularly when it comes to higher throughput needs, increasing the time and costs of research.

SignalStar mIHC technology has established an assay that circumvents these issues, providing a sensitive, reproducible, and customizable assay for comprehensive biomarker spatial imaging of fixed tissues. Obtaining results is 70% faster than other mIHC approaches with no need for time consuming assay optimization, and results are obtained in two days from start to finish, speeding up translational workflows.

In addition to being a trusted leader of high quality and rigorously tested antibodies, CST provides comprehensive technical support and educational assistance available for the set-up and follow through of your imaging experiments to ensure their successful completion and streamlining your research process.



For Research Use Only. Not for Use in Diagnostic Procedures.