

Highly Multiplexed IHC Assays to Examine Immune Checkpoints and Biomarkers for Immunotherapy

INTRODUCTION

The emergence of an increasing number of immunotherapy biomarkers and the importance of their context within the tumor microenvironment has resulted in a need for high-plex immunohistochemistry (IHC) assays. Using highly specific and validated antibodies developed for this purpose, we constructed several fluorescent multiplexed, TSA-based assays to examine the frequency, spatial localization, and proximity of immune cells within the tumor microenvironment.

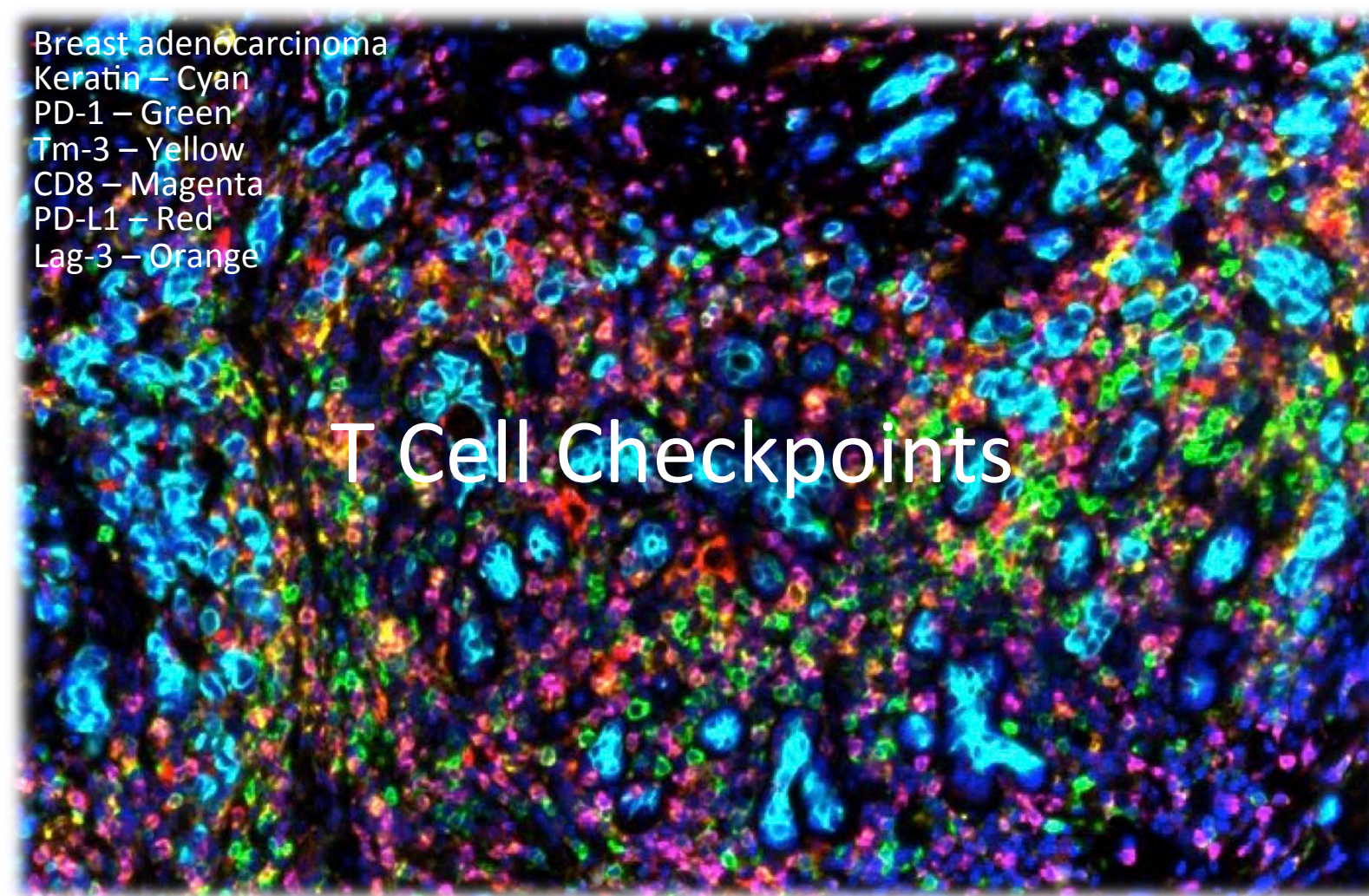
Our data demonstrates the feasibility of simultaneous detection of seven fluorochromes in order to visualize immunosuppressive receptors associated with the exhausted T cell phenotype, myeloid-derived suppressor cells, and the PD-1:PD-L1 axis. Our findings demonstrate the utility of multiplex IHC to deconvolute protein expression and interactions within the complex tumor microenvironment.

METHODS

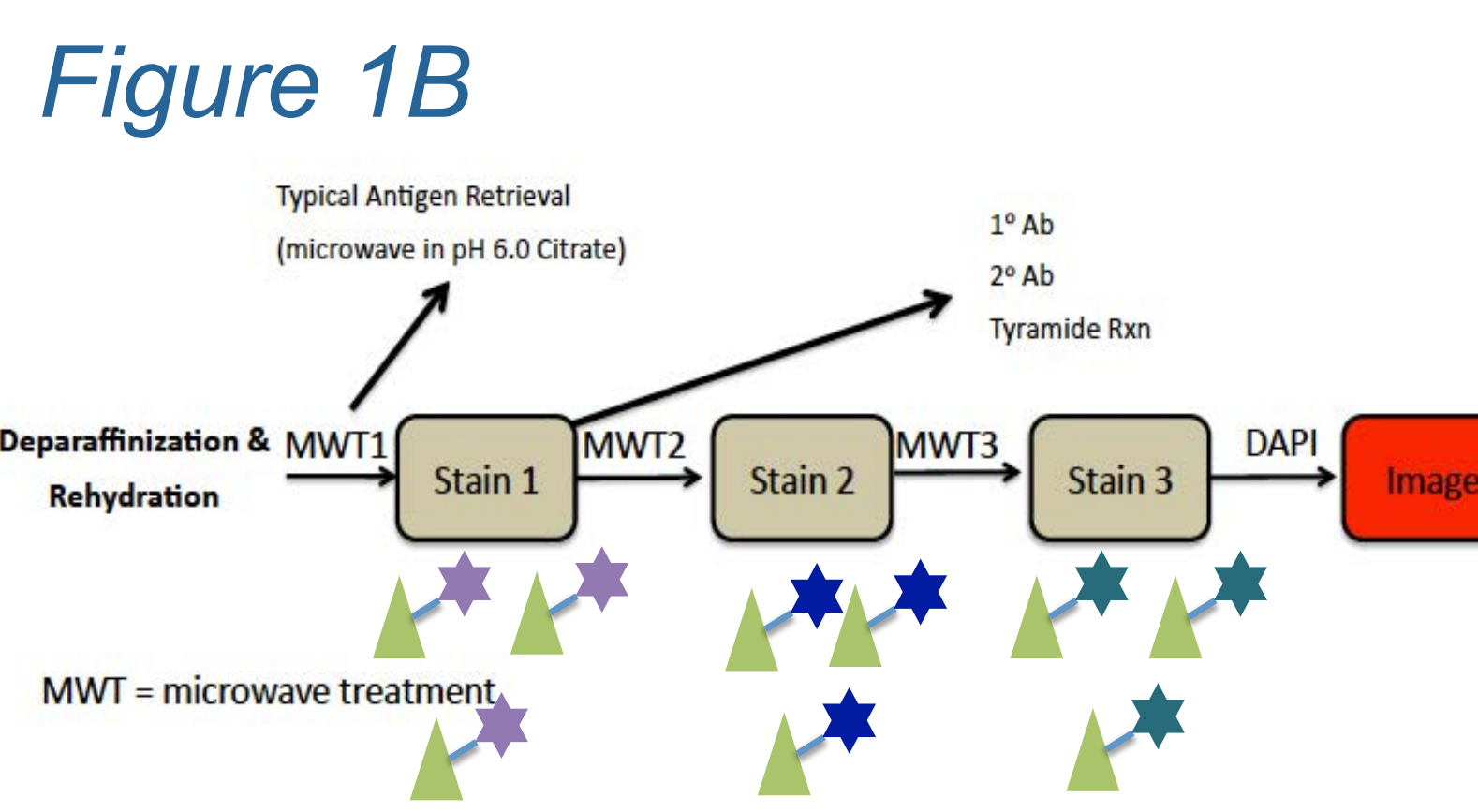
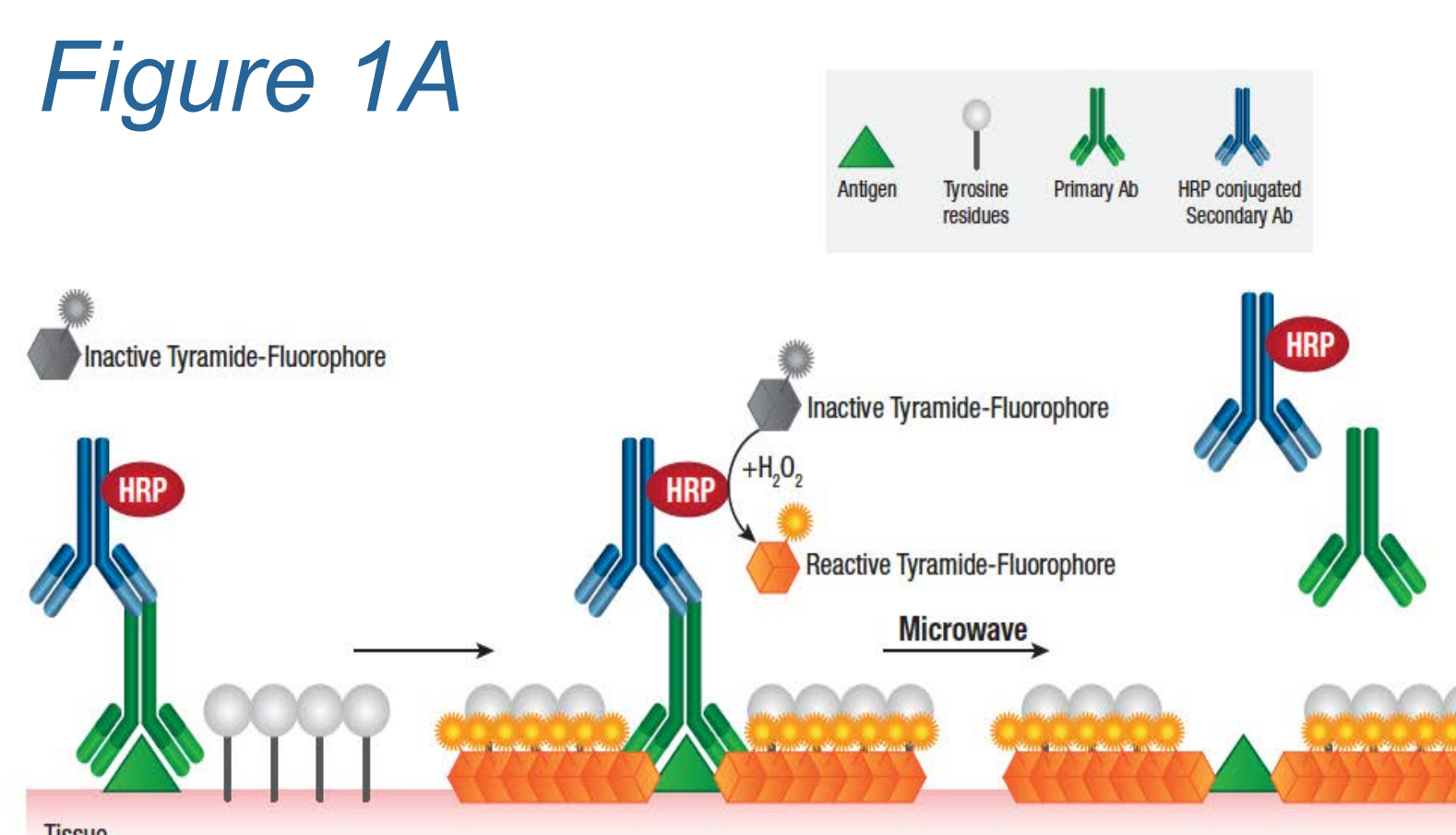
Tyramide signaling amplification (TSA) was used to serial stain tumor tissue of various types. This protocol allows for the use of multiple rabbit monoclonal antibodies in a single panel. A Mantra quantitative pathology workstation (PerkinElmer) was used to spectrally unmix the fluorescent signal in each image, and the InForm Image Analysis software (PerkinElmer) was used to provide quantitative data. Immuno-oncology-centric panels have been constructed, as well as those that focus on receptors involved in certain targeted cancer therapies.

CONCLUSIONS

- Multiplex IHC panels consisting of up to six targets plus DAPI were constructed and validated in various tumor types.
- Highly detailed images illustrating the utility of mIHC to detect:
 - Spatial localization of immune cells within the tumor micro-environment.
 - Co-localization and frequency of immune checkpoint receptors.
 - Proximity of suppressive immune cells and their interactions with tumor cells in the process of immune evasion.
- Any Cell Signaling Technology, Inc. IHC-validated antibody can be used to construct mIHC panels.



Seven-plex IHC of key immunology phenotypic markers and therapeutic targets



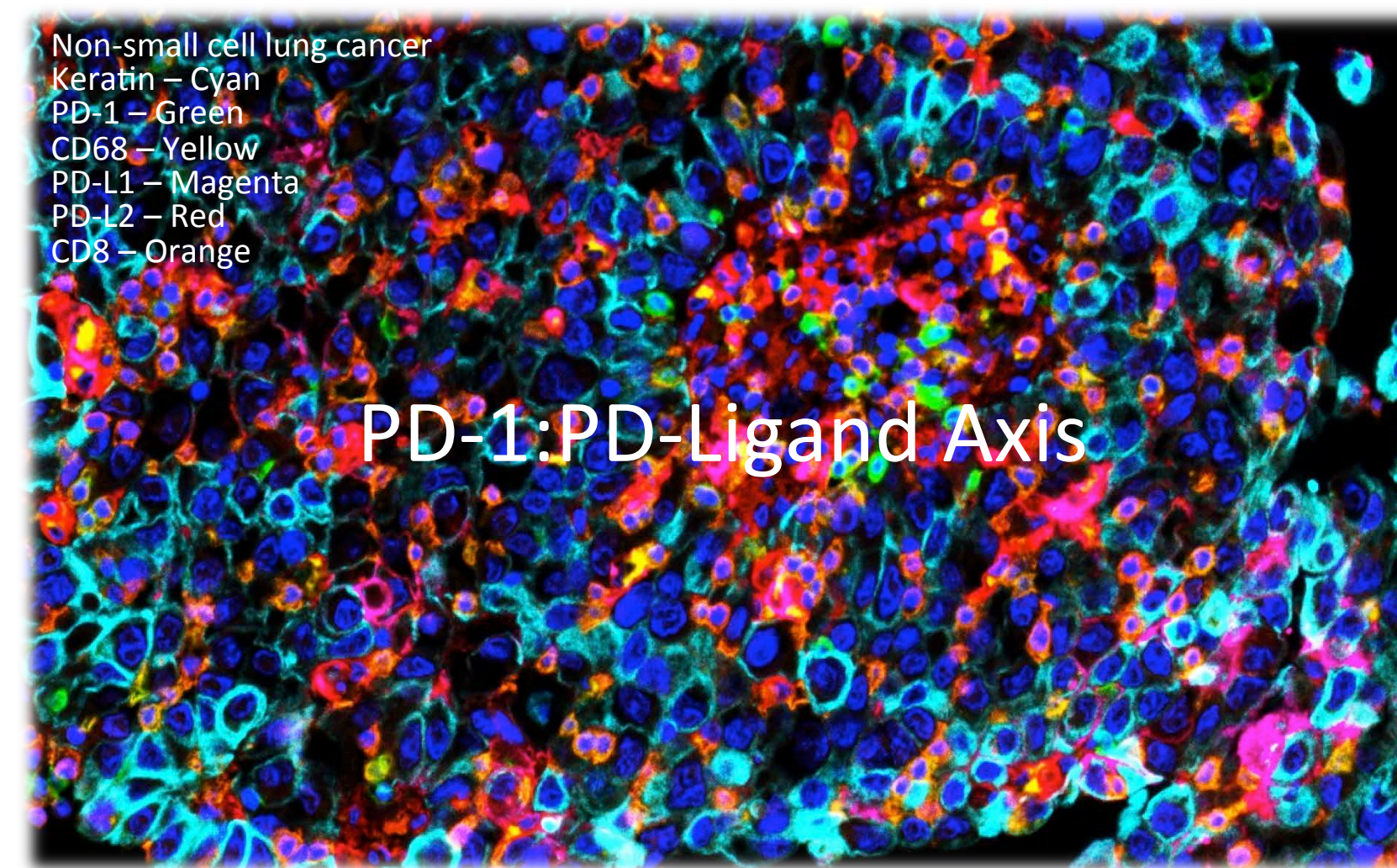
Targeted Therapies	Panel	Targets
Xalkori	ALK (D5F3B) XP® Rabbit mAb #3633 ROS1 (D4D6B) Rabbit mAb #15027 Met (D1C2) XP® Rabbit mAb #8198 Krt (C11) Mouse mAb #4545	ALK (D5F3B) XP® Rabbit mAb #3633
		ROS1 (D4D6B) Rabbit mAb #15027
		Met (D1C2) XP® Rabbit mAb #8198
Prostate	PSA (D11E1) XP® Rabbit mAb #2475 PSMA (D7I8E) XP® Rabbit mAb #12815 AR (D6F1) XP® Rabbit mAb #5153 Krt (C11) Mouse mAb #4545	PSA (D11E1) XP® Rabbit mAb #2475
		PSMA (D7I8E) XP® Rabbit mAb #12815
		AR (D6F1) XP® Rabbit mAb #5153
EGFR	EGFR (D38B1) XP® Rabbit mAb #4267 Her-2 (D8F12) XP® Rabbit mAb #4290 Her-3 (D22C5) XP® #12708 Krt (C11) Mouse mAb #4545	EGFR (D38B1) XP® Rabbit mAb #4267
		Her-2 (D8F12) XP® Rabbit mAb #4290
		Her-3 (D22C5) XP® #12708

Immunotherapy	Panel	Targets
Effector Cell/T _H Ratio	CD8a (C8/144B) Mouse mAb #70306 FoxP3 (D2W8E™) Rabbit mAb #98377 Krt (C11) Mouse mAb #4545	CD8a (C8/144B) Mouse mAb #70306
		FoxP3 (D2W8E™) Rabbit mAb #98377
		Krt (C11) Mouse mAb #4545
Co-regulatory Ligand/T-cell Panel	PD-L1 (E1L3N®) XP® Rabbit mAb #13684 B7-H4 (D1M8) XP® Rabbit mAb #14572 CD8 (C8/144B) Mouse mAb #70306 FoxP3 (D2W8E™) Rabbit mAb #98377 Krt (C11) Mouse mAb #4545	PD-L1 (E1L3N®) XP® Rabbit mAb #13684
		B7-H4 (D1M8) XP® Rabbit mAb #14572
		CD8 (C8/144B) Mouse mAb #70306
Her-2/Co-Regulatory Ligand Co-expression	PD-L1 (E1L3N®) XP® Rabbit mAb #13684 B7-H3 (D9M2L) XP® #14058 B7-H4 (D1M8) XP® Rabbit mAb #14572 Her-2 (D8F12) XP® Rabbit mAb #4290 Krt (C11) Mouse mAb #4545	PD-L1 (E1L3N®) XP® Rabbit mAb #13684
		B7-H3 (D9M2L) XP® #14058
		B7-H4 (D1M8) XP® Rabbit mAb #14572
Anti-Mouse	PD-L1 (E1L3N®) XP® Rabbit mAb #13684 CD-8 (C8/144B) Mouse mAb #70306 Tim-3 (D5D5R™) XP® Rabbit mAb #45208 Arginase-1 (D4E3M™) XP® #93688 Krt (C11) Mouse mAb #4545	PD-L1 (E1L3N®) XP® Rabbit mAb #13684
		CD-8 (C8/144B) Mouse mAb #70306
		Tim-3 (D5D5R™) XP® Rabbit mAb #45208
Exhausted T-cell	Tim-3 (D5D5R™) XP® Rabbit mAb #45208 PD-1 (D4W2J) XP® Rabbit mAb #86163 CD8 (C8/144B) Mouse mAb #70306 Lag-3 (D2G4D™) XP® Rabbit mAb #15372 CD68 (D4B9C) XP® #76437 Krt (C11) Mouse mAb #4545	Tim-3 (D5D5R™) XP® Rabbit mAb #45208
		PD-1 (D4W2J) XP® Rabbit mAb #86163
		CD8 (C8/144B) Mouse mAb #70306
PD-1:PD-L1 Ligand Axis	PD-L1 (E1L3N®) XP® Rabbit mAb #13684 PD-1 (D4W2J) XP® Rabbit mAb #86163 CD68 (D4B9C) XP® #76437 CD8 (C8/144B) Mouse mAb #70306 Tim-3 (D5D5R™) XP® Rabbit mAb #45208 Krt (C11) Mouse mAb #4545	PD-L1 (E1L3N®) XP® Rabbit mAb #13684
		PD-1 (D4W2J) XP® Rabbit mAb #86163
		CD68 (D4B9C) XP® #76437

Figure 1A. A schematic of the Tyramide Signaling Amplification (TSA) protocol. FFPE tissue is deparaffinized and rehydrated. After antigen retrieval, tissue is stained with a 1^o antibody followed by an HRP conjugated 2^o antibody. The HRP enzyme catalyzes the reaction of tyramide (PerkinElmer) to a reactive form, which binds tyrosine residues on and near the target. The tyramide is conjugated to a fluorophore which can then be detected and imaged.
Products used:
SignalStain® Boost IHC Detection Reagent (HRP, Rabbit) #8114
SignalStain® Boost IHC Detection Reagent (HRP, Mouse) #8125
SignalStain® Antibody Diluent #8112
SignalStain® Mounting Medium #14177

Figure 1B. A schematic illustrating the TSA protocol as applied to the serial staining process necessary for multiplexing. To multiplex, the tissue is then heated such that the weaker hydrogen bonds formed by the primary and secondary antibodies are broken, but the covalent bonds between the deposited tyramide and tyrosine residues remain. Multiplex staining is achieved in a serial fashion, with another staining cycle started once the first primary/secondary antibody pair has been removed.

Table 1A. A list of the mIHC panels that have been constructed to date, grouped according to their research area. CST's highly specific and validated antibodies serve as the foundation for clarifying distinct signals.



Spatial localization and frequency of exhausted T cells expressing immune checkpoint receptors

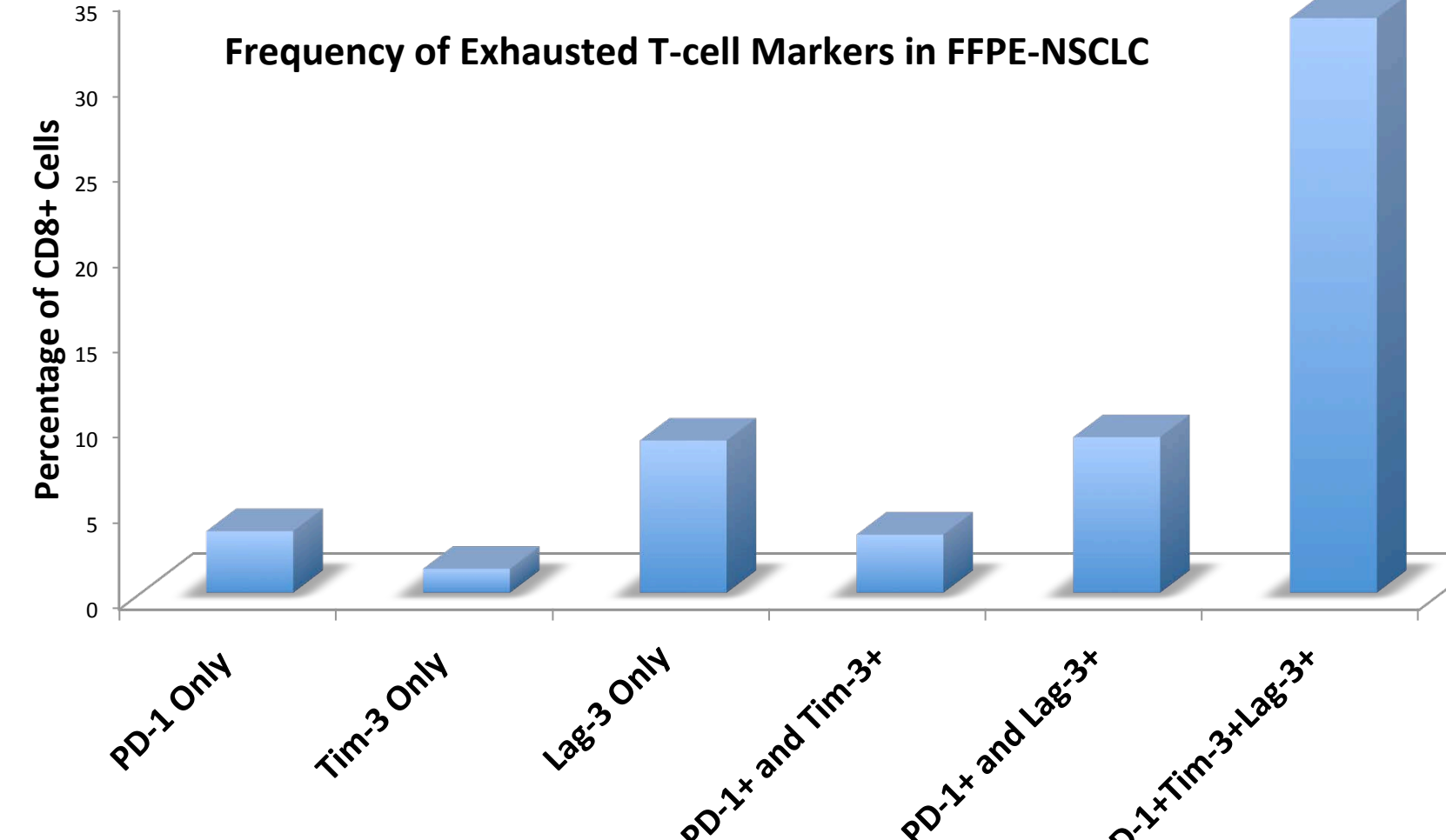
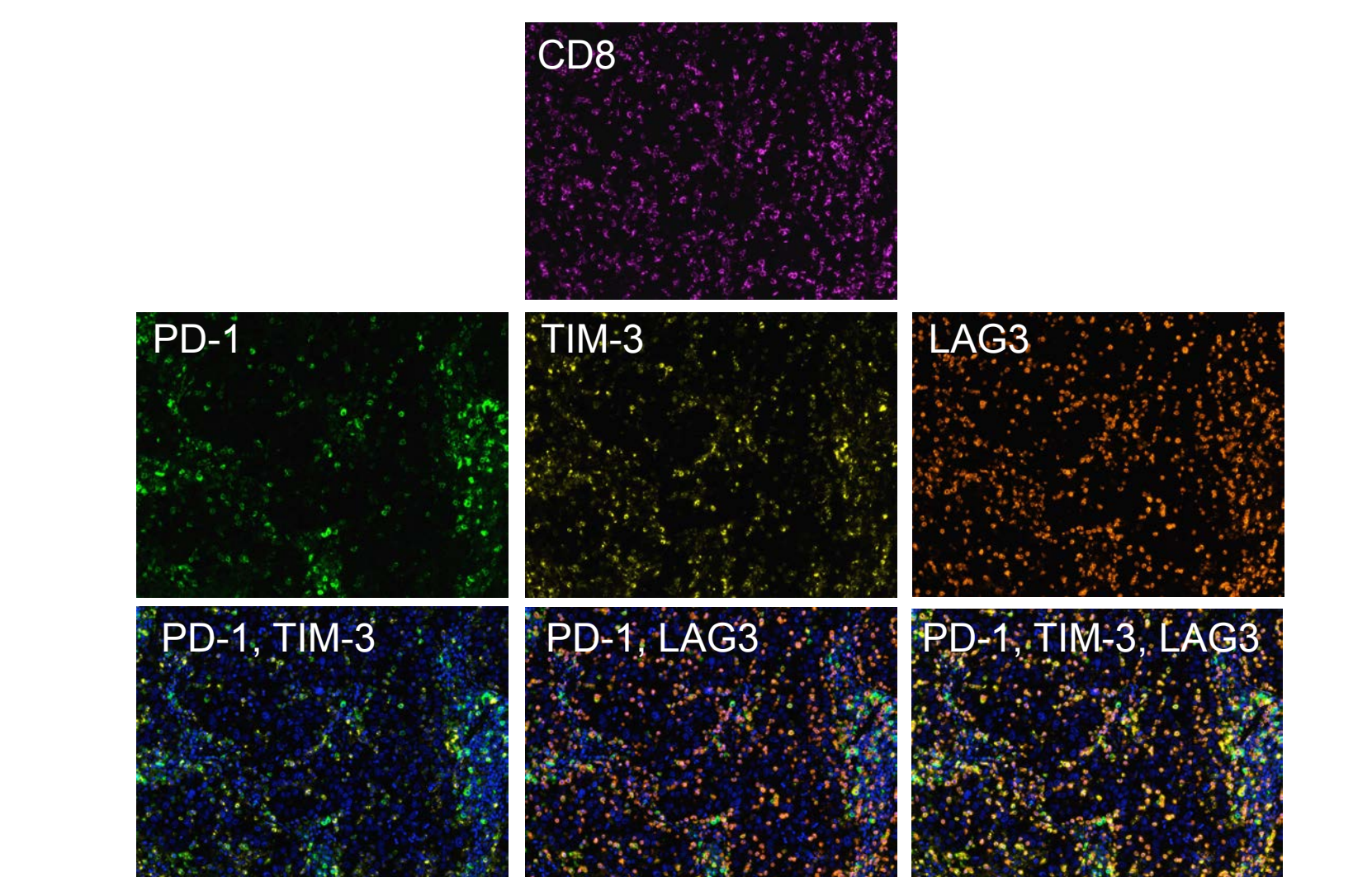
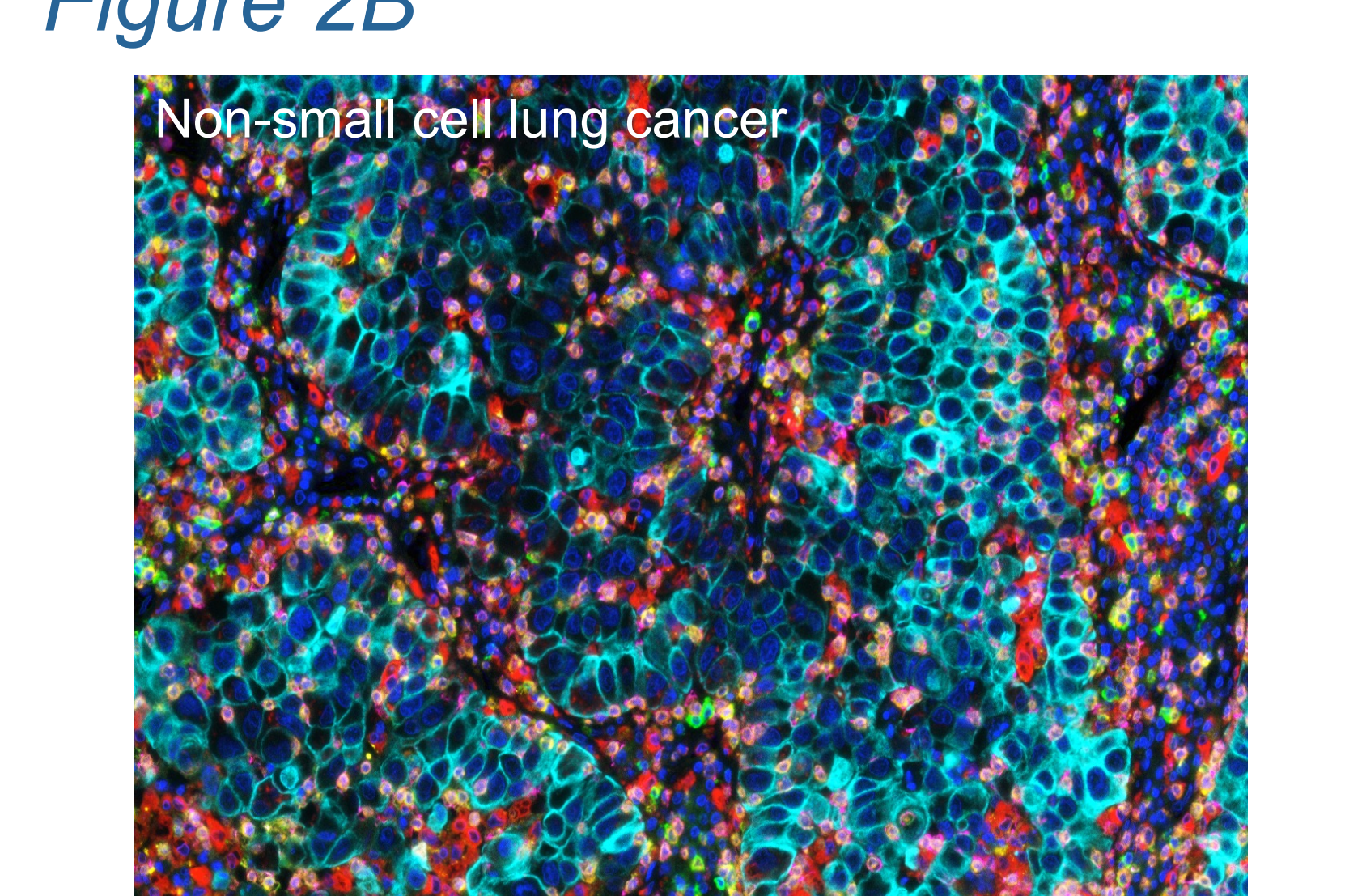
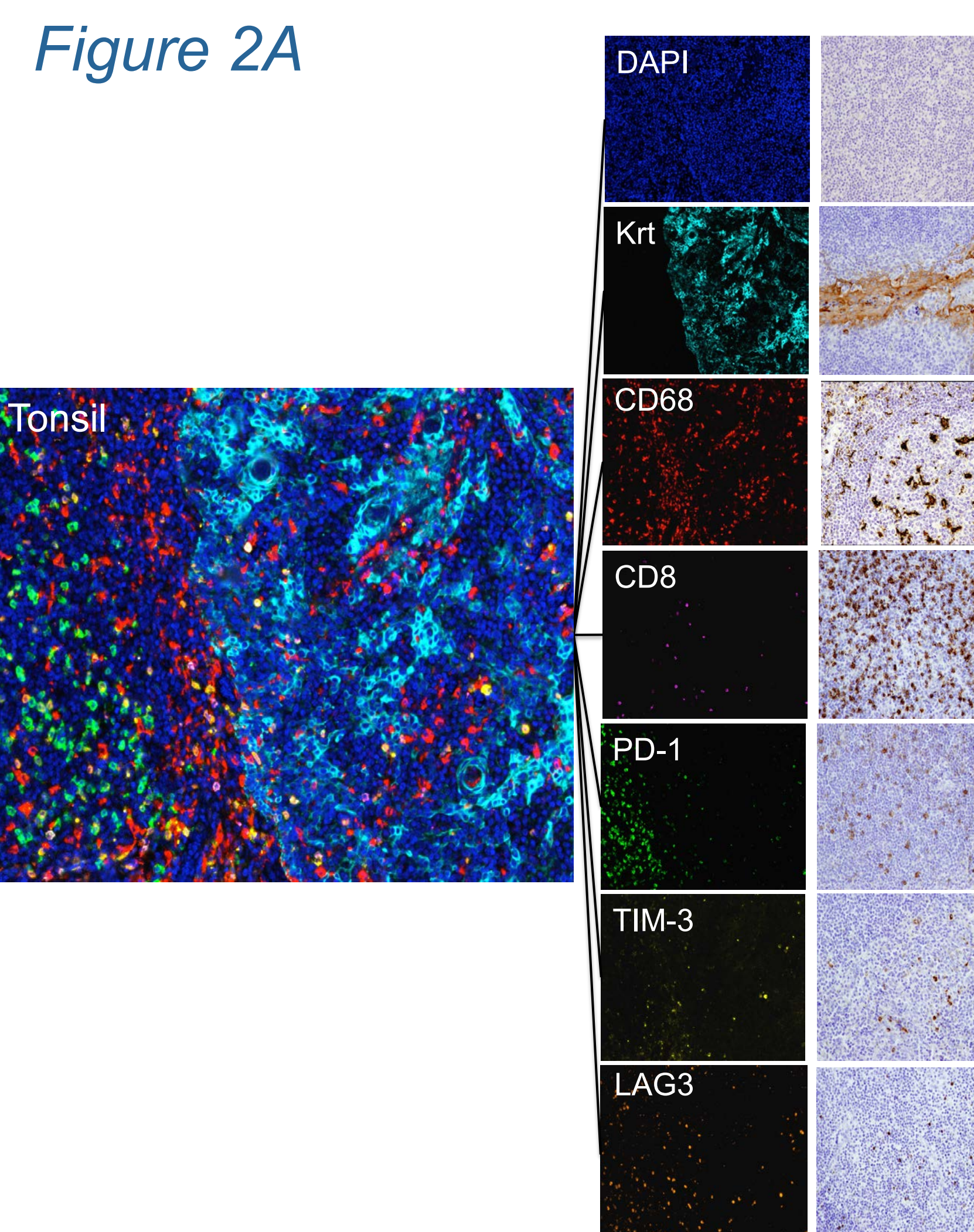
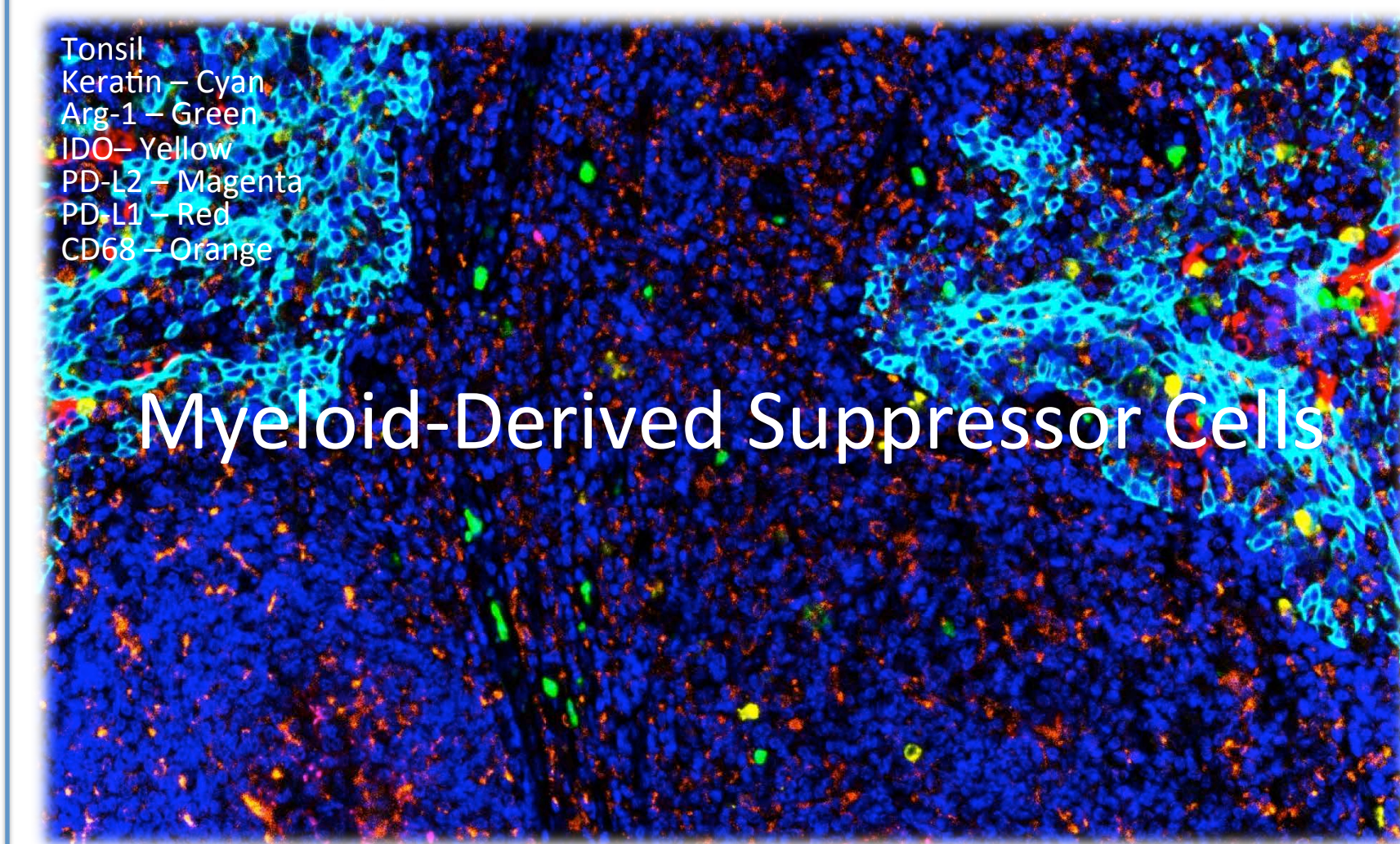


Figure 2A. Normal human tonsil was stained with an exhausted T cell phenotype mIHC panel demonstrating seven fluorescent signals can be clearly unmixed with spectral imaging. The fluorescent multiplex staining in each case mirrors the corresponding chromogenic staining performed on serial sections from the same tissue block.

Figure 2C. A NSCLC section was stained with an exhausted T cell phenotype mIHC panel including PD-1, TIM-3 and LAG3. Co-expression of these markers was imaged and quantified.



Proximity and co-localization of immunosuppressive cells within the tumor microenvironment

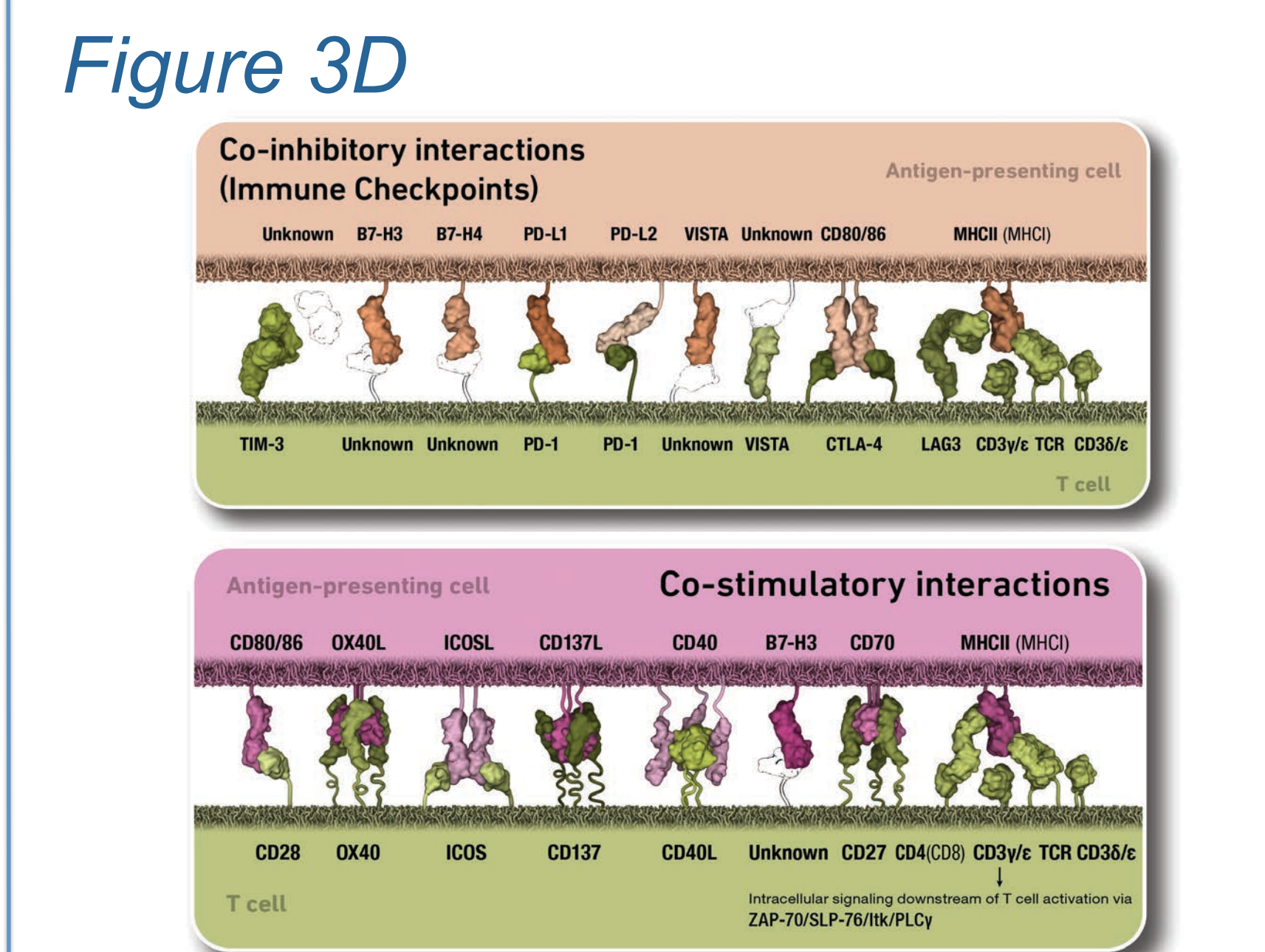
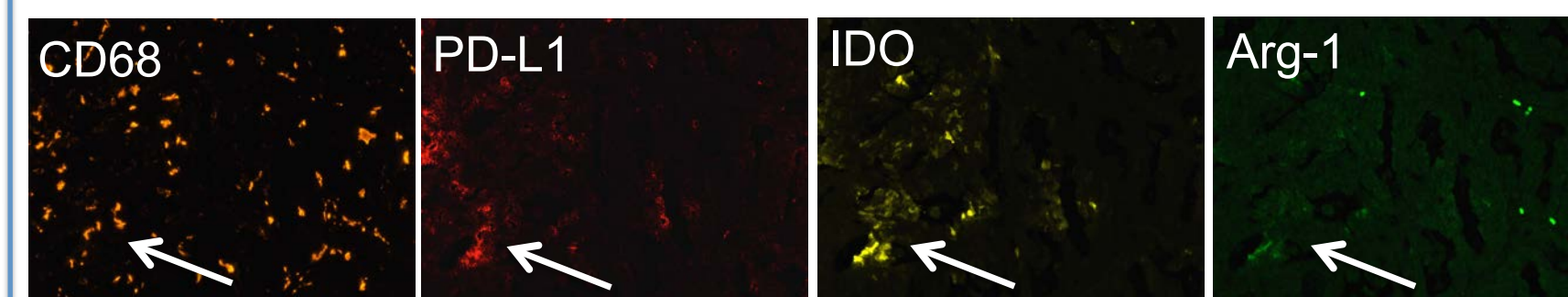
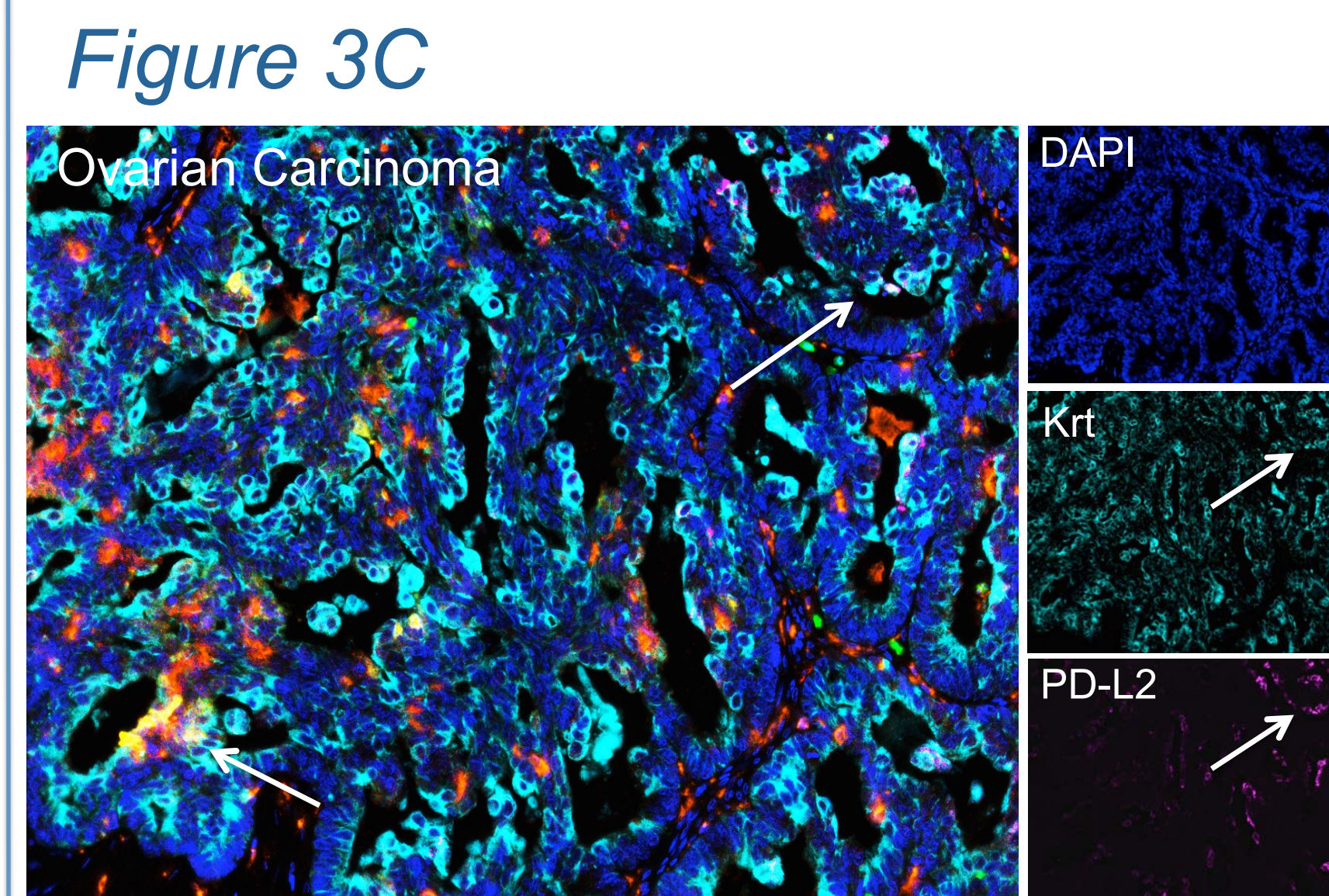
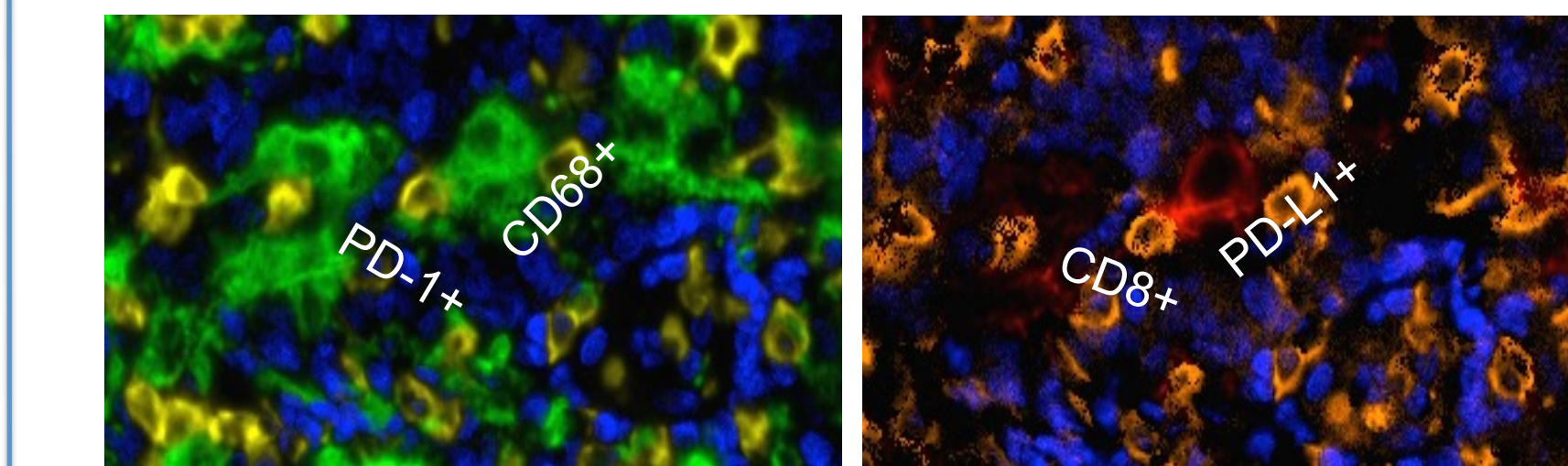
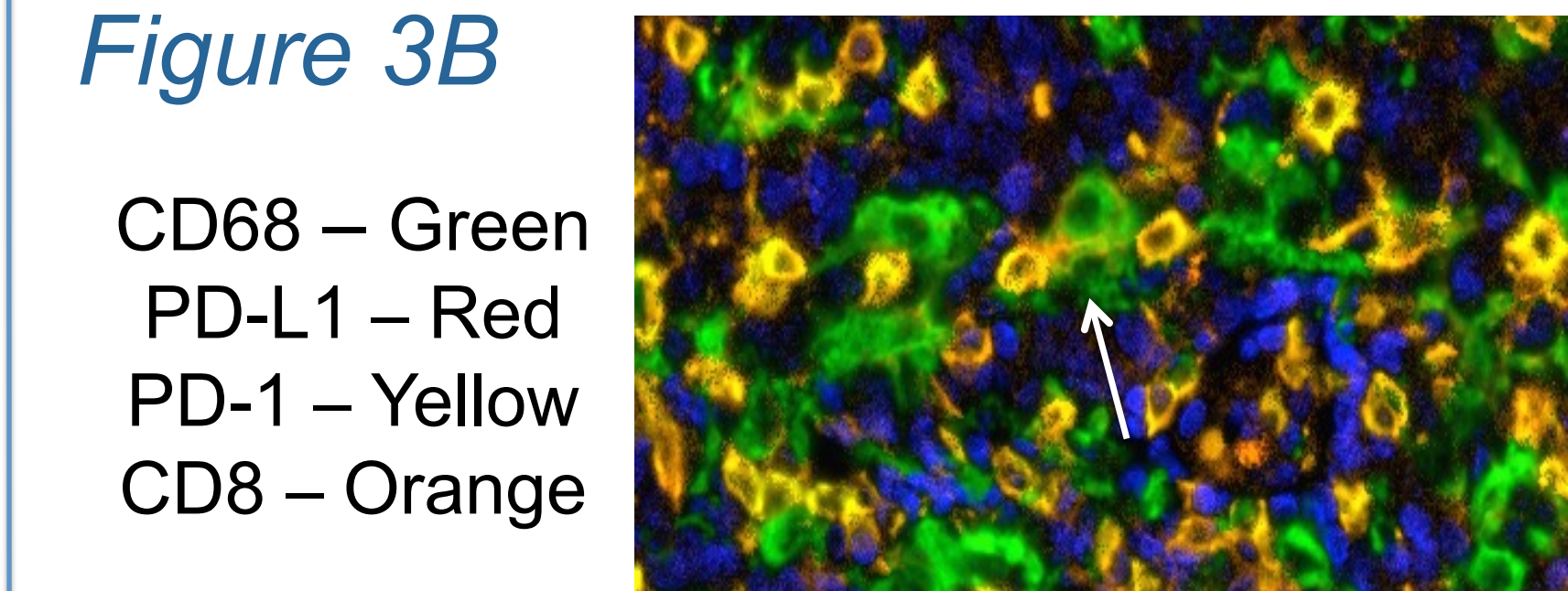
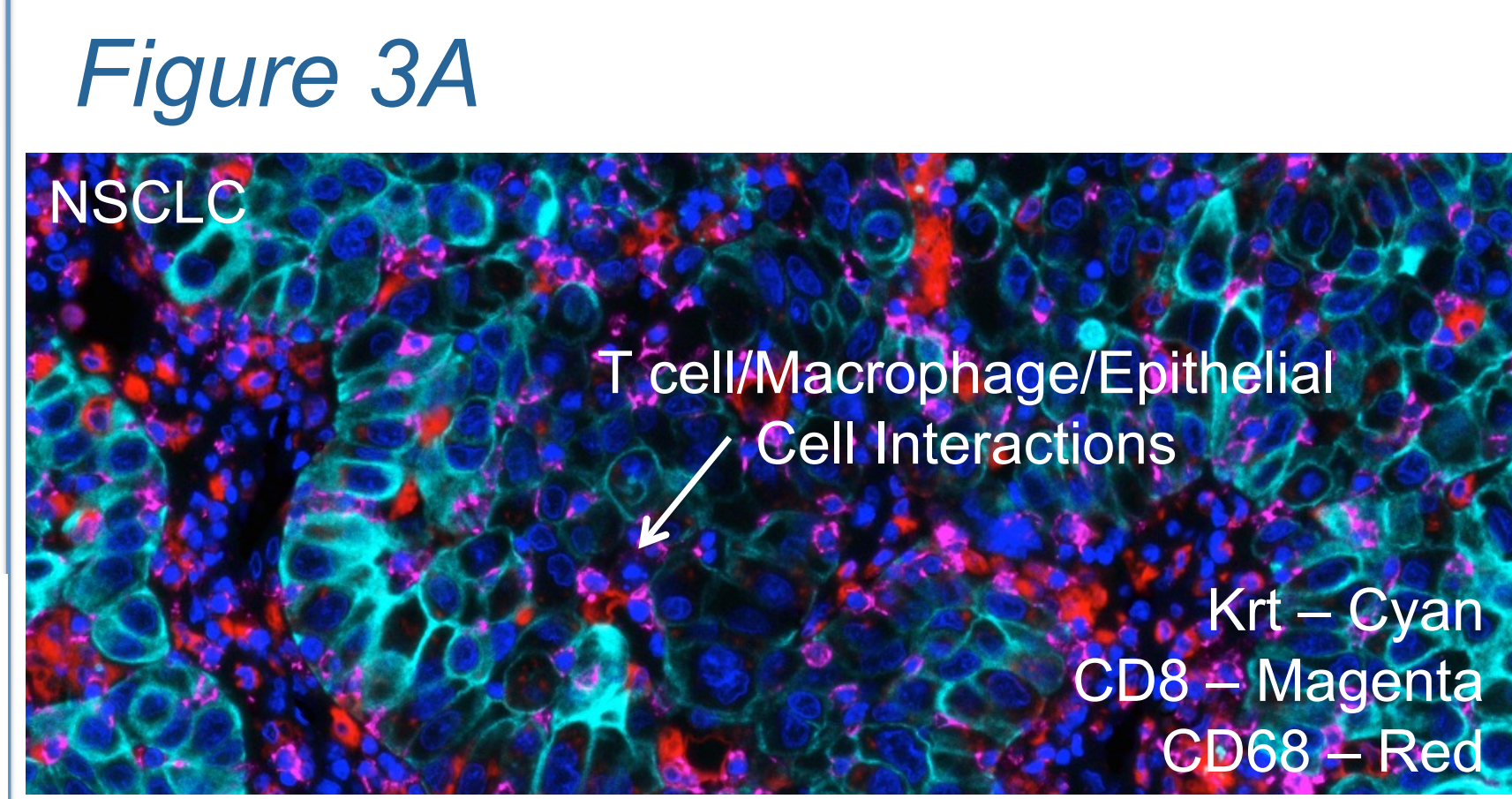


Figure 3A. CD8+ T cells and CD68+ macrophages are shown in close proximity to tumor cells in a NSCLC section, suggesting possible interactions.

Figure 3B. PD-L1+CD68+ macrophages interacts with a PD-1+CD8+ T cells in FFPE B cell Non-Hodgkin's lymphoma.

Figure 3C. Ovarian carcinoma was stained with a MDCS-centric mIHC panel. Co-expression of PD-L1, IDO, and Arginase-1 was detected in CD68+ macrophages. PD-L2 co-localized with Krt+ tumor cells.

Figure 3D. Future panels will continue to incorporate various co-inhibitory and co-stimulatory receptors that have been identified as being important in tumor immune evasion.