

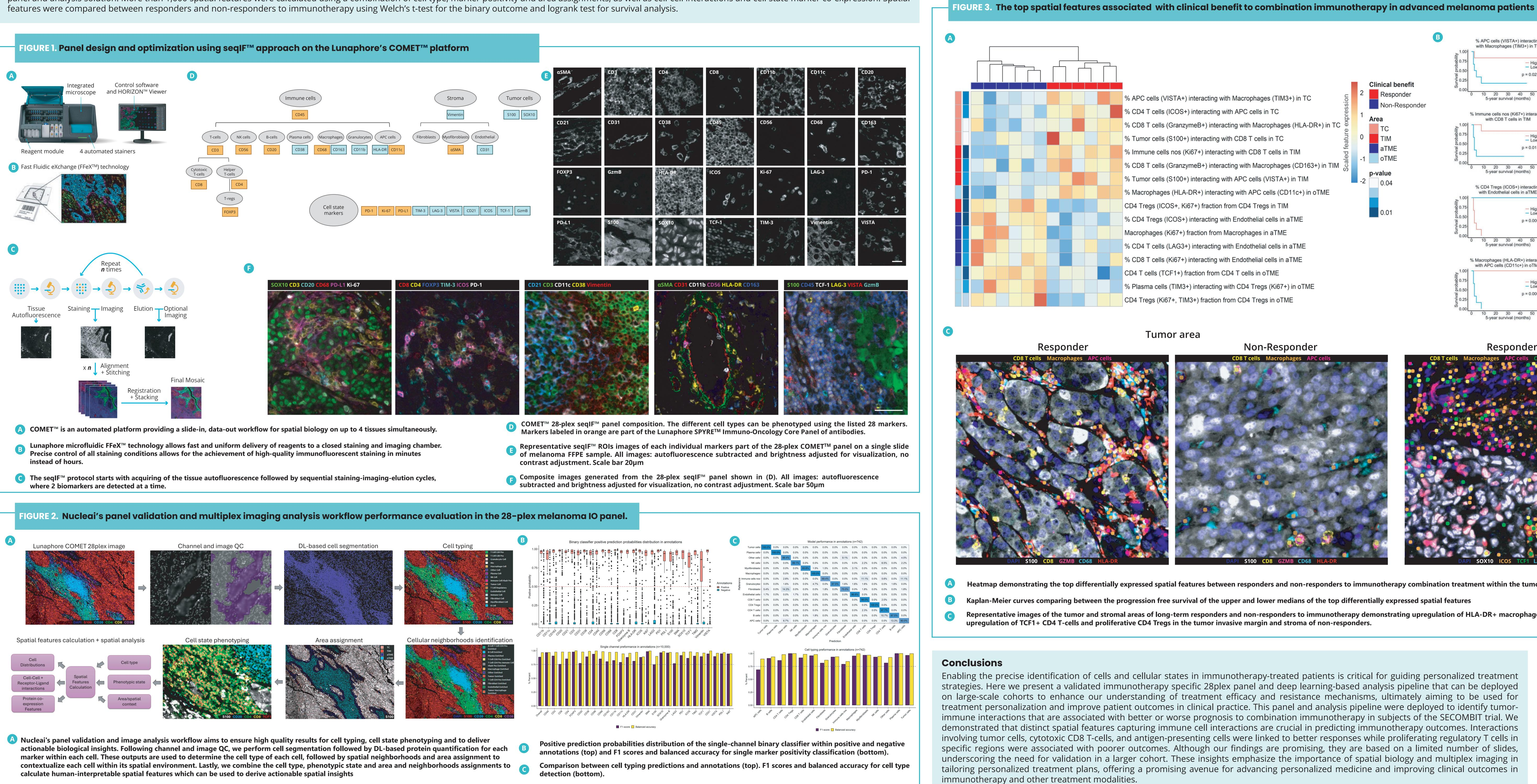
# Application of a Novel Multiplex Imaging-based Immunotherapy Panel and Al-powered Analysis Solution for Spatial Biomarker Identification on Immunotherapy-treated Melanoma Patients

Ettai Markovits<sup>2</sup>, Marion Bonnet<sup>1</sup>, Pedro Machado Almeida<sup>1</sup>, Gabriele Madonna<sup>3</sup>, Mariaelena Capone<sup>3</sup>, Michael Mints<sup>2</sup>, Paolo Ascierto<sup>3,4</sup>, Antonio Sorrentino<sup>1</sup> Sorrentino<sup>1</sup>, Saska Brajkovic<sup>1</sup>, Oscar Puig<sup>2</sup>, Paolo Ascierto<sup>3,4</sup>, Antonio Sorrentino<sup>1</sup> <sup>1</sup>Lunaphore, Tolochenaz, Switzerland<sup>2</sup>Nucleai, Tel Aviv, Israel<sup>3</sup>Istituto Nazionale Tumori IRCCS Fondazione "G. Pascale Napoli", Napoli, Italy<sup>4</sup>On behalf of SECOMBIT study group<sup>4</sup>

## Introduction

Identifying biomarkers that predict patient response to immunotherapy is critical for optimizing treatment strategies and improving clinical outcomes. Despite the success of immunotherapy, a significant proportion of patients do not respond to treatment. Thus, there is an urgent need for more robust methods to differentiate responders. Spatial biology, which examines the precise localization and interaction of cells within the tissue, plays a crucial role in understanding the complex tumor-immune microenvironment. In this study, we present a novel multiplex imaging-based immunotherapy panel and a comprehensive analysis pipeline to characterize the spatial distribution and function of immune cells and its application in a cohort of immunotherapy- and targeted therapy-treated melanoma patients.

### Methods



## We designed a 28-plex panel to perform sequential immunofluorescence (seqIF<sup>M</sup>) on the COMET<sup>M</sup> platform (Figure 1A-C) [1] to target key biomarkers associated with tumor microenvironment composition (TME), immune cell infiltration, and immune checkpoint pathways (Figure 1D-F). Utilizing Nucleai's deep-learning-based multiplex imaging analysis pipeline (Figure 2A) [2], we identified 16 cell types, including 10 different immune cell populations, in addition to 10 cell state markers. Single channel and cell typing, cellular neighborhoods were designated as previously described [3], and were assigned to the tumor or stromal areas (Figure 2B-C). The tumor area was further divided into the tumor core (TC) and tumor invasive margin (TIM), while the stroma was divided into adjacent TME (aTME) and outer TME (oTME) based on the distance from the tumor-stroma border. We obtained pre-treatment biopsies from patients in arm B (Ipilimumab + Nivolumab until progression, followed by BRAF targeted therapy (Encorafenib + Binimetinib)) from the SECOMBIT Phase II Trial (NCT02631447) [4, 5] with known long-term response or rapid progression to immunotherapy combination treatment. These samples were profiled using the aforementioned panel and analysis solution. More than 1,000 spatial features were calculated using a combination of cell type, marker positivity and area assignments, as well as cell-cell interactions and cell state marker co-expression. Spatial

## Results

Our novel multiplex imaging panel and analysis pipeline demonstrated high accuracy in both cell typing and quantification of protein expression as demonstrated by a balanced accuracy (> 0.9) in all markers and F1 scores (> 0.75) in >97% of markers, as well as F1 scores > 0.75 in >90% of cell types. Combination of cell types. Combination of cell types cell phenotypic state and the spatial context of the cells enables the quantification of known biomarkers such as T cell activation states, T cell infiltration patterns, and TLS maturation, as well as discovery of novel biomarkers that capture biologically relevant cell states and tumor-immune cells and immune-immune cells interactions. Comparison of calculated spatial features between long-term responders (n=6) and rapid progressors (progressors (progressors (progressors (progressors (progressors (progressors)) and rapid progressors (progressors) and activation status across the tumor areas associated with immunotherapy response. Within the tumor cells, cytotoxic CD8 T-cells and antigen presenting cells (APC) such as HLA-DR+ macrophages and dendritic cells were associated with better immunotherapy outcome. In contrast, a high percentage of proliferating regulatory T cells within the tumor invasive margin (TIM) was associated with a worse clinical outcome from combination immunotherapy. In the adjacent TME (aTME), endothelial cell interactions with T-cells and macrophage proliferation were associated with immunotherapy resistance, while in the outer TME (oTME), regulatory Tcell proliferation as well as regulatory T-cells – plasma cell interactions and TCF-1 positivity within CD4 T-cells were associated with lack of response. In contrast, the interaction between MHC class II macrophages and APC cells was associated with improved clinical outcome.

Taken together, we demonstrate that area-specific immunotherapy response and highlight the importance of spatial biology in predicting immunotherapy outcomes and tailoring personalized treatment strategies.

Heatmap demonstrating the top differentially expressed spatial features between responders and non-responders to immunotherapy combination treatment within the tumor core (TC), tumor invasive margin (TIM), adjacent TME (aTME) and outer TME (oTME) Kaplan-Meier curves comparing between the progression free survival of the upper and lower medians of the top differentially expressed spatial features Representative images of the tumor and stromal areas of long-term responders and non-responders and APC, and Granzyme B+ CD8 T-cells in the tumor area of responders and

Enabling the precise identification of cells and cellular states in immunotherapy-treated patients is critical for guiding personalized treatment strategies. Here we present a validated immunotherapy specific 28plex panel and deep learning-based analysis pipeline that can be deployed on large-scale cohorts to enhance our understanding of treatment efficacy and resistance mechanisms, ultimately aiming to be used for treatment personalization and improve patient outcomes in clinical practice. This panel and analysis pipeline were deployed to identify tumorimmune interactions that are associated with better or worse prognosis to combination immunotherapy in subjects of the SECOMBIT trial. We demonstrated that distinct spatial features capturing immune cell interactions are crucial in predicting immunotherapy outcomes. Interactions involving tumor cells, cytotoxic CD8 T-cells, and antigen-presenting cells were linked to better responses while proliferating regulatory T cells in specific regions were associated with poorer outcomes. Although our findings are promising, they are based on a limited number of slides, underscoring the need for validation in a larger cohort. These insights emphasize the importance of spatial biology and multiplex imaging in tailoring personalized treatment plans, offering a promising avenue for advancing personalized medicine and improving clinical outcomes in

Scan to get a digital version of the poster

Tumore core % APC cells (VISTA+) interacting % CD4 T cells (ICOS+) interacting % CD8 T cells (GranzymeB+) interacting % Tumor cells (S100+) interacting with Macrophages (TIM3+) in TC with APC cells in TC with Macrophages (HLA-DR+) in TO with CD8 T cells in TC High
Low High
 Low p = 0.0270 = 0.0210 20 30 40 50 5-year survival (months) 5-year survival (months) 5-year survival (months 5-year survival (months) Non-Responder Tumor invasive margin % CD8 T cells (GranzymeB+) interacting % Tumor cells (S100+) interacting CD4 Tregs (ICOS+, Ki67+) fraction from % Immune cells nos (Ki67+) interacting with CD8 T cells in TIM with APC cells (VISTA+) in TIM CD4 Tregs in TIM with Macrophages (CD163+) in TIM - High 10 20 30 40 50 5-year survival (months) 20 30 40 50 5-year survival (months) 5-year survival (months) 5-year survival (months) % CD4 T cells (LAG3+) interactir % CD8 T cells (Ki67+) interacting 6 CD4 Tregs (ICOS+) interacting with Endothelial cells in aTME with Endothelial cells in aTME with Endothelial cells in aTME Macrophages in aTME p = 0.0004910 20 30 40 50 10 20 30 40 50 20 30 40 5-year survival (months) 5-year survival (months) 5-year survival (months) 5-year survival (months) **Outer TME** CD4 T cells (TCF1+) fraction from % Macrophages (HLA-DR+) interacting CD4 Treas (Ki67+, TIM3+) fraction fro % Plasma cells (TIM3+) interacting with APC cells (CD11c+) in oTME CD4 T cells in oTME with CD4 Tregs (Ki67+) in oTME CD4 Tregs in oTME - High - Low p = 0.000490 = 0.0004910 20 30 40 50 20 30 40 50 20 30 40 50 60 5-year survival (months 5-vear survival (months 5-vear survival (months

**Tumor Invasive Margin and Stroma** 

Responder

# References

- . Rivest F, et al. Fully automated sequential immunofluorescence (seqIF) for hyperplex spatial proteomics. Sci Rep. 2023, 13(1):16994.
- Markovits E, et al. A novel deep learning pipeline for cell typing and phenotypic marker
- quantification in multiplex imaging, bioRxiv 2022;
- S. Schürch, C. M. et al. Coordinated Cellular Neighborhoods Orchestrate Antitumoral Immunity at the Colorectal Cancer Invasive Front. Cell 182, 1341-1359.e19 (2020).
- 4. Ascierto PA, et al. Sequencing of Ipilimumab Plus Nivolumab and Encorafenib Plus Binimetinib for Untreated BRAF-Mutated Metastatic Melanoma (SECOMBIT): A Randomized, Three-Arm, Open-Label Phase II Trial. J Clin Oncol. 2023 Jan 10;41(2):212-221
- Ascierto PA, et al. Sequential immunotherapy and targeted therapy for metastatic BRAF V600 mutated melanoma: 4-year survival and biomarkers evaluation from the phase II SECOMBIT trial. Nat Commun. 2024, 15(1):146
- . Ascierto PA, et al. Sequencing of checkpoint of BRAF/MEK inhibitors on brain metastasis in Melanoma. NEJM Evid 2024;3(10)

