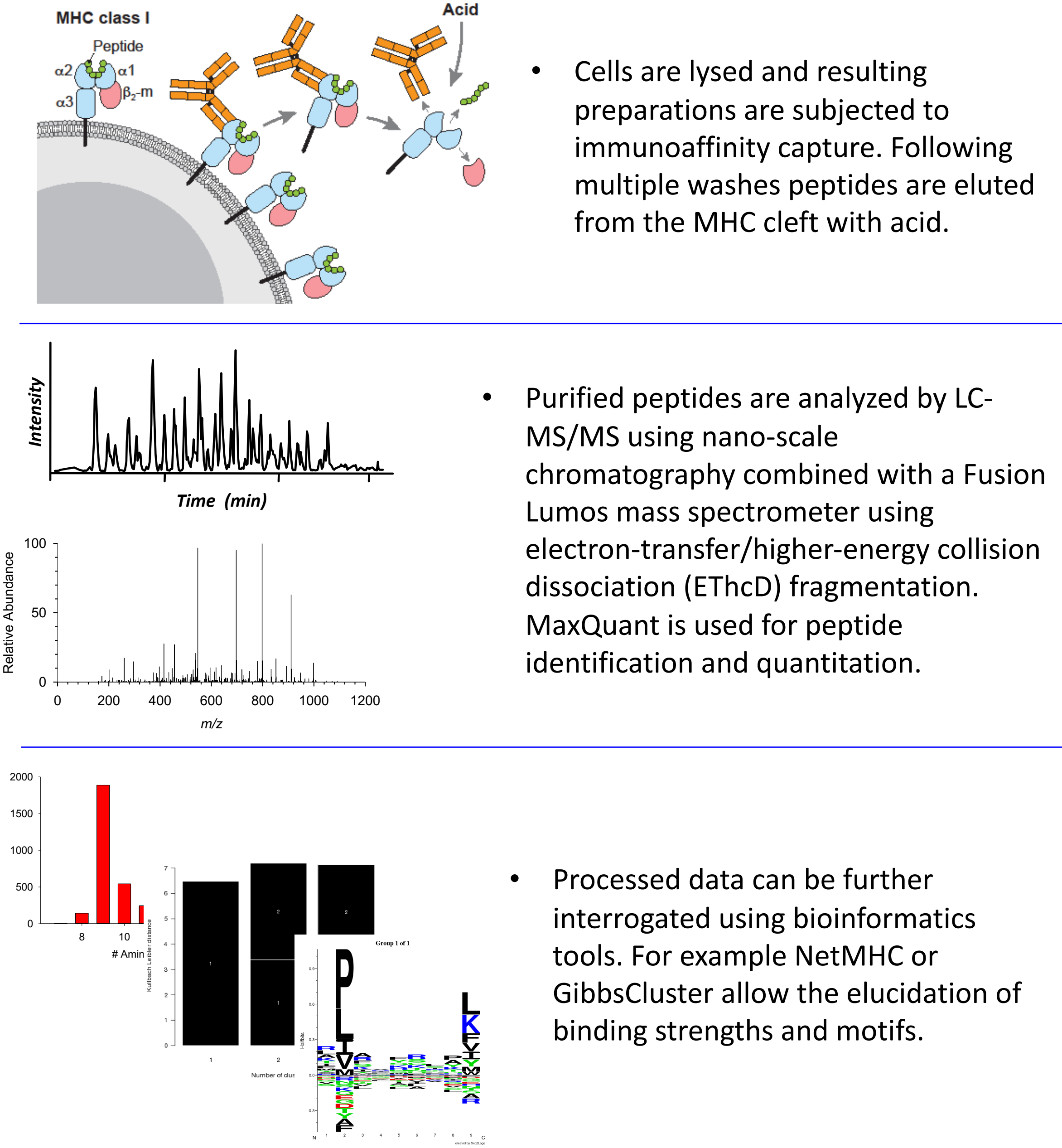
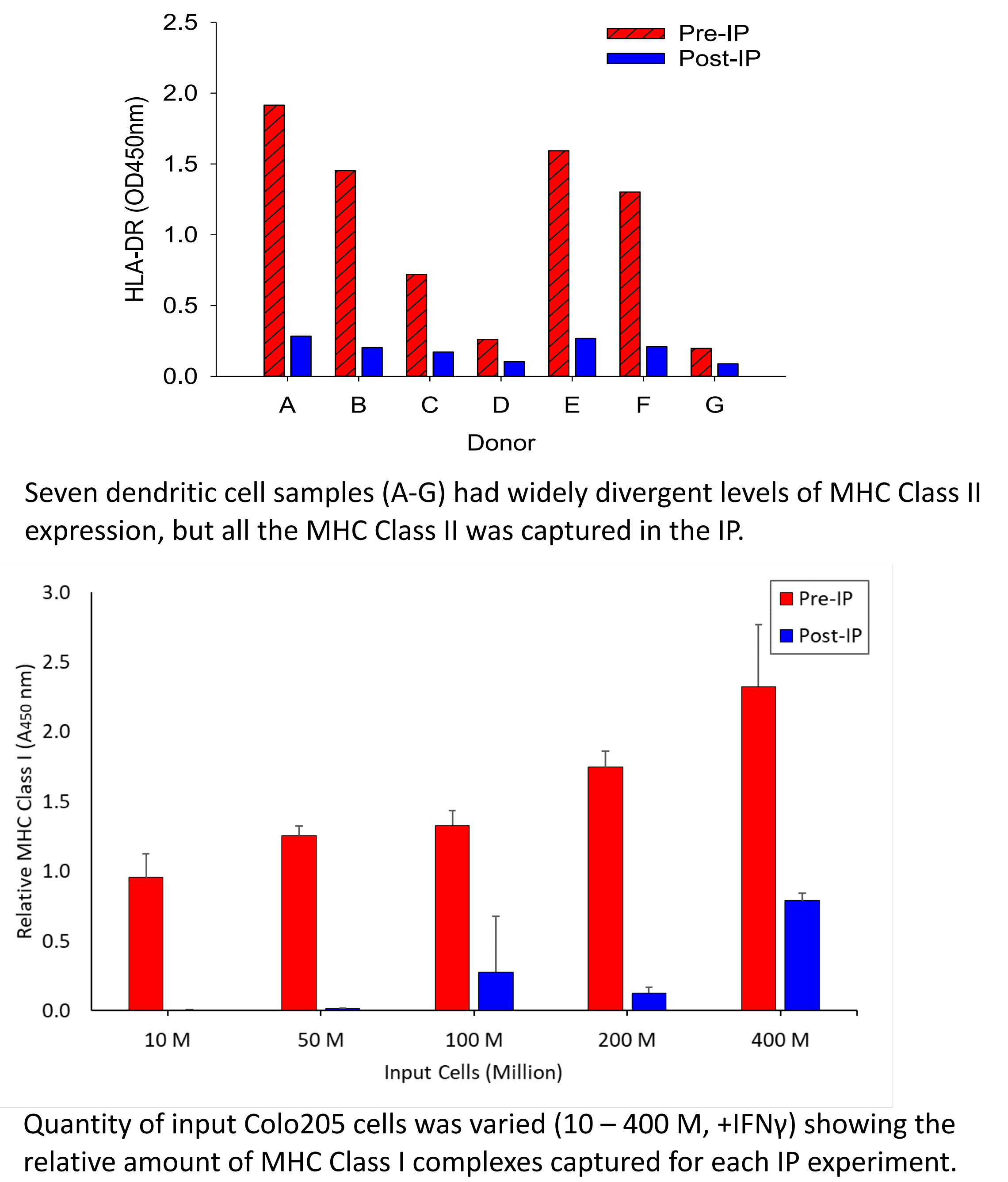


Workflow

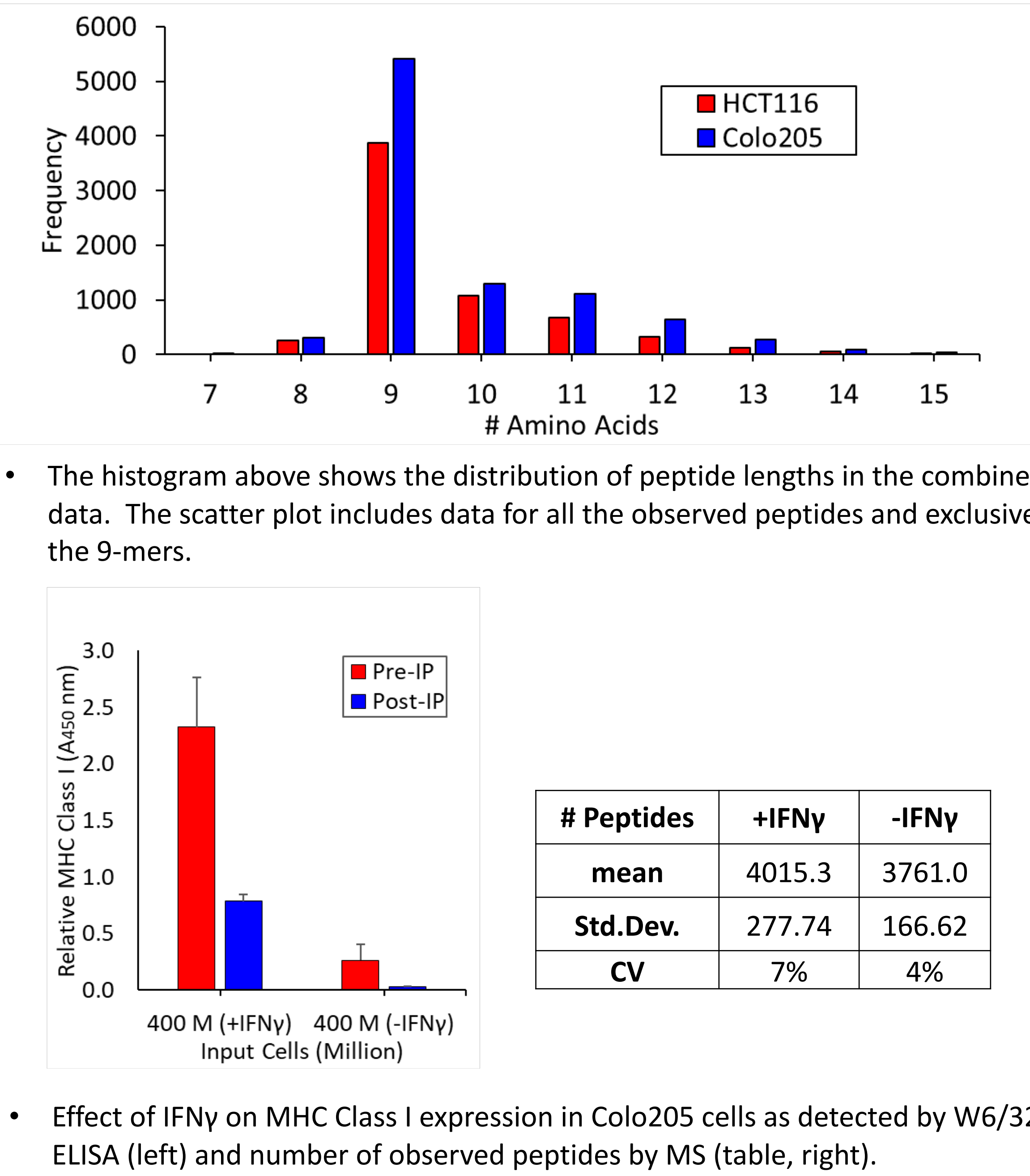
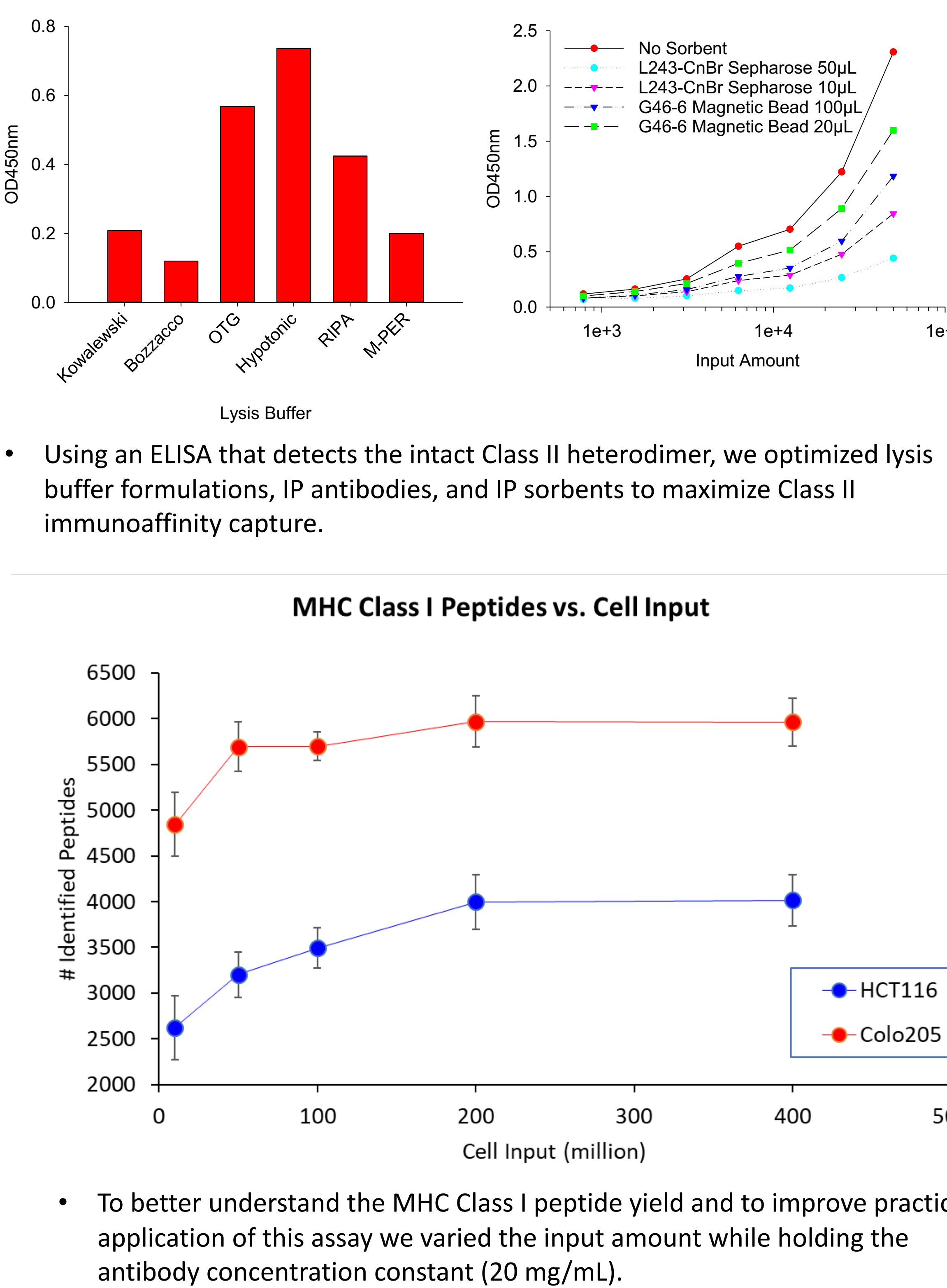


MHC Recovery

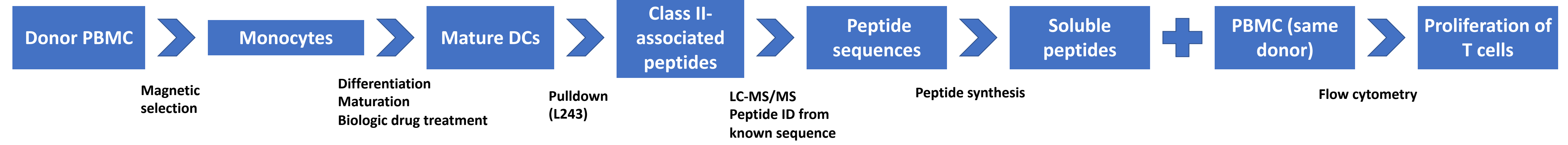
- Custom ELISAs detect intact MHC Class I and MHC Class II complexes to monitor recovery of the MHC molecules.
- This is used for both tissue/cell specific optimization and as a quality control check during the studies.



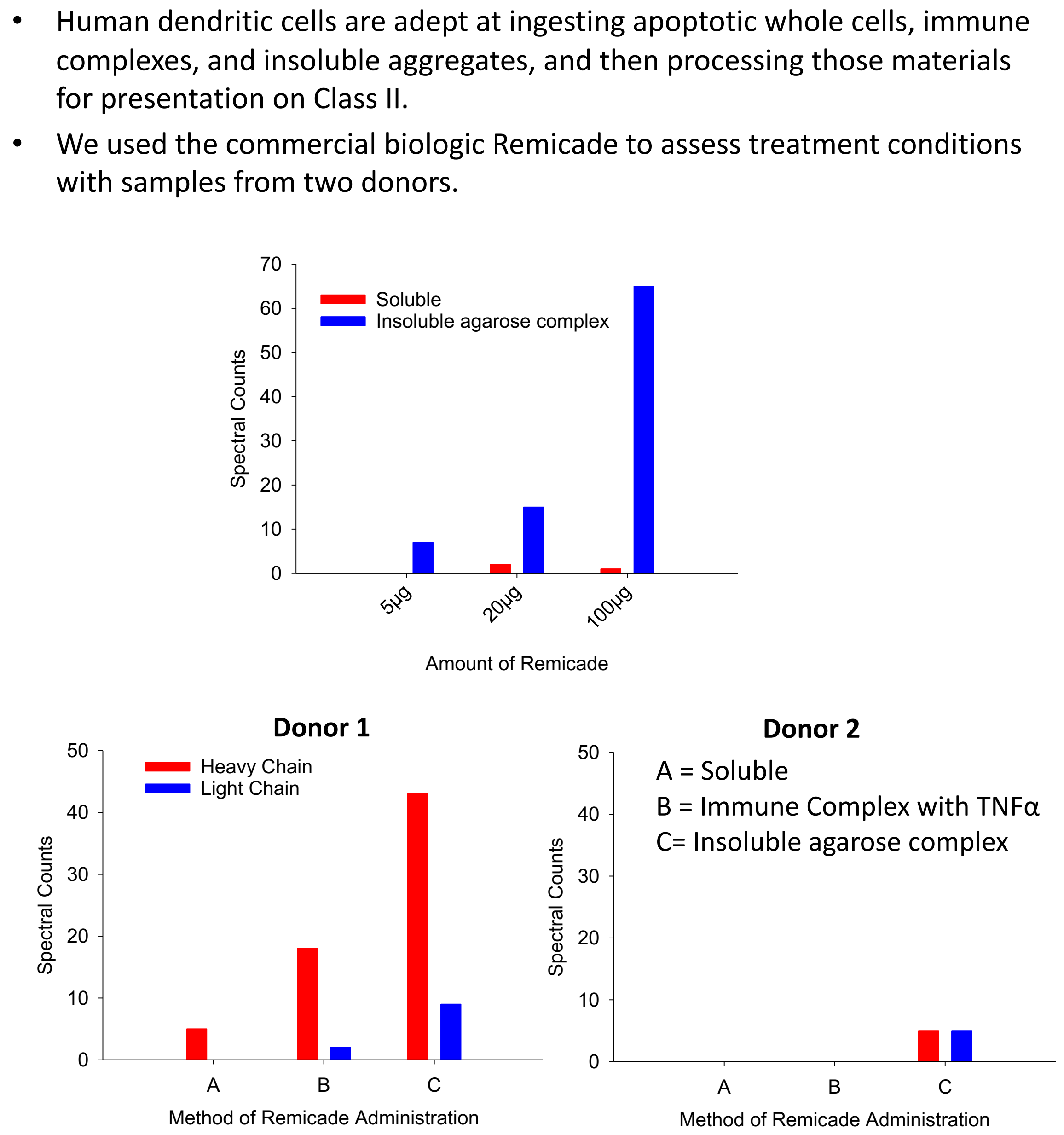
Optimization



Biotherapeutic T-Cell Epitopes

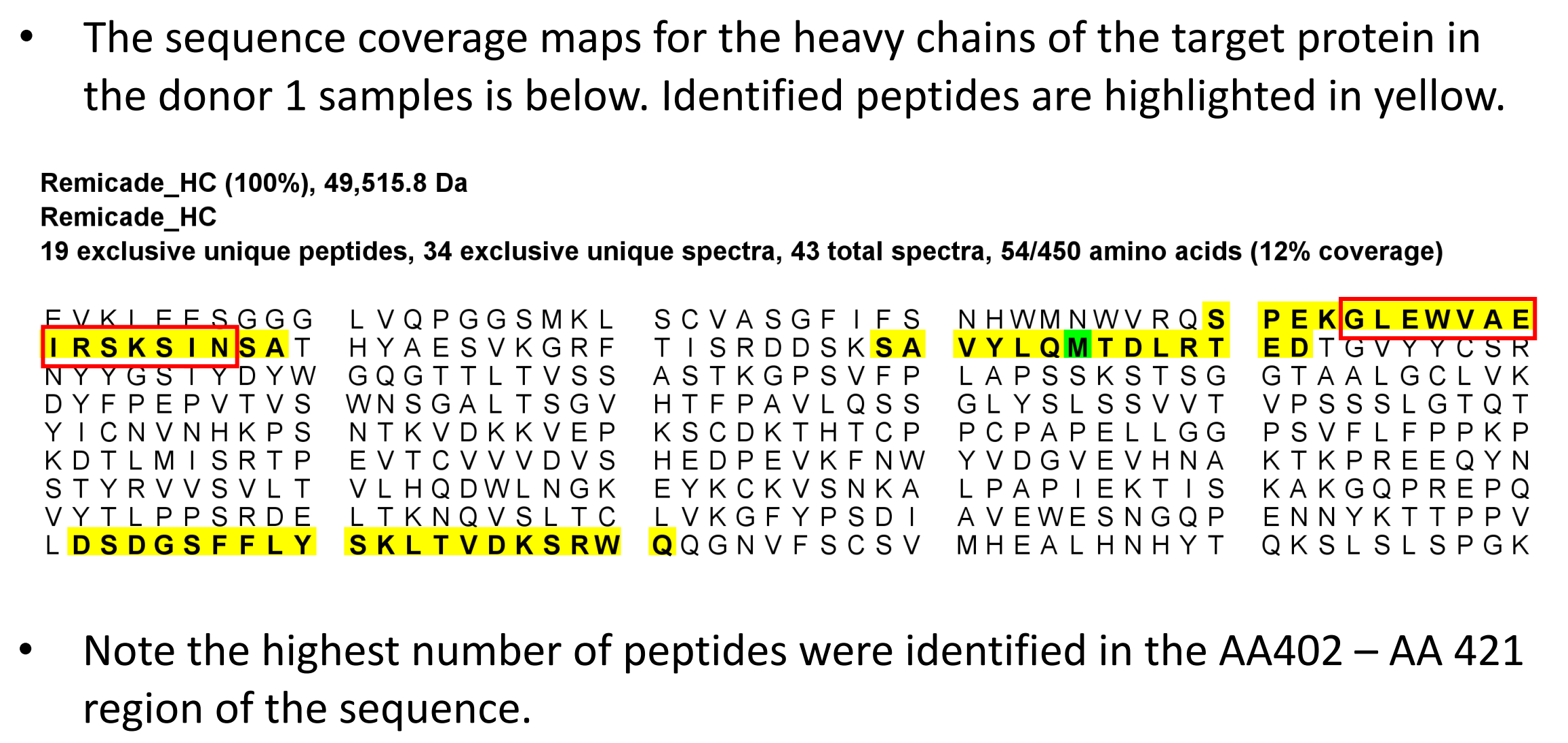


DC Pulsing With Biotherapeutic



- Dendritic cells prefer immune complexes or insoluble aggregates over soluble proteins for ingestion, processing, and display on Class II.
- Soluble protein antigen (Remicade) is taken up and processed far less efficiently than aggregated antigen.
- The ability to present peptides derived from Remicade is donor-specific. Increasing the concentration of soluble protein antigen does little to improve the outcome.

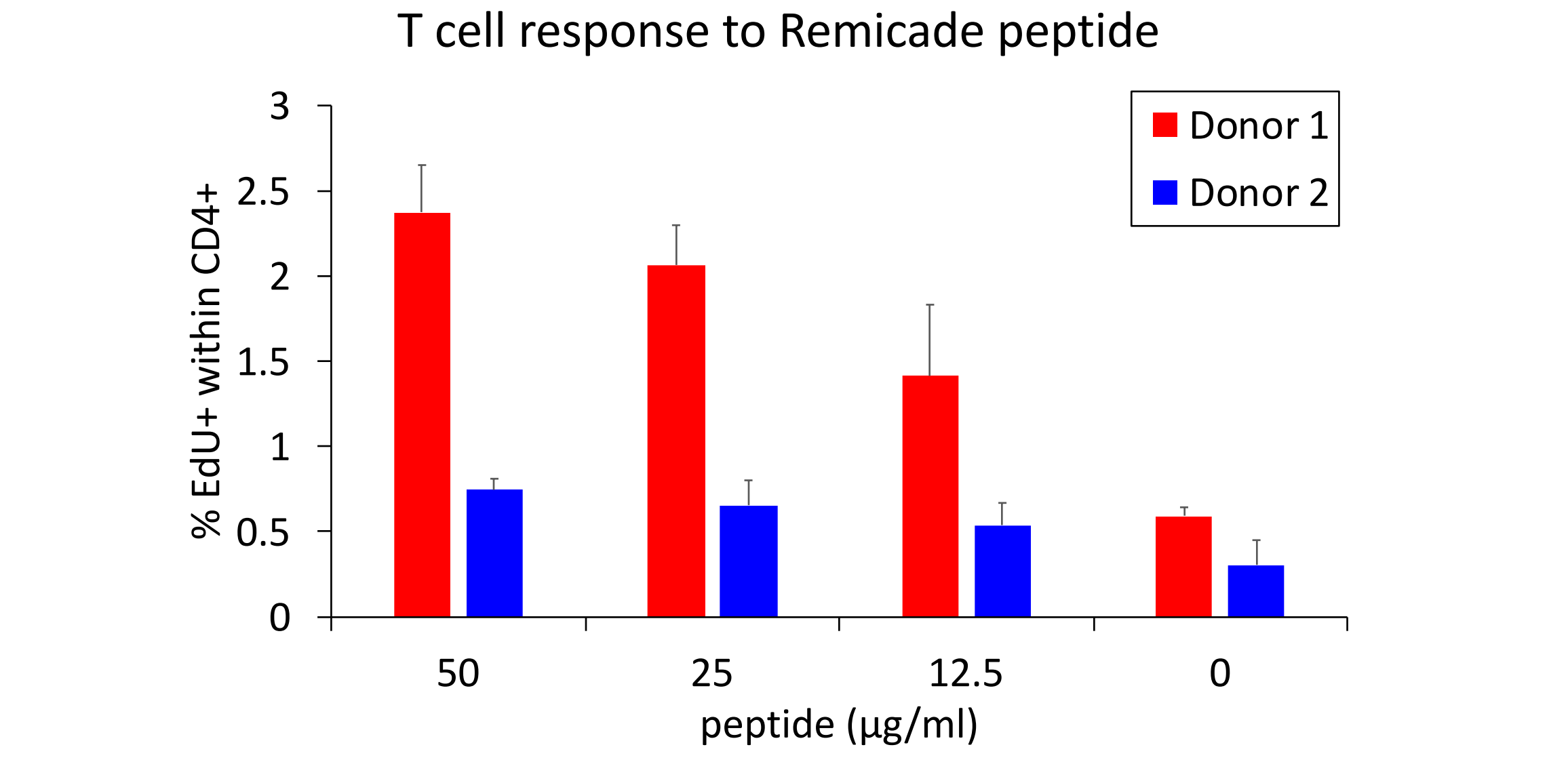
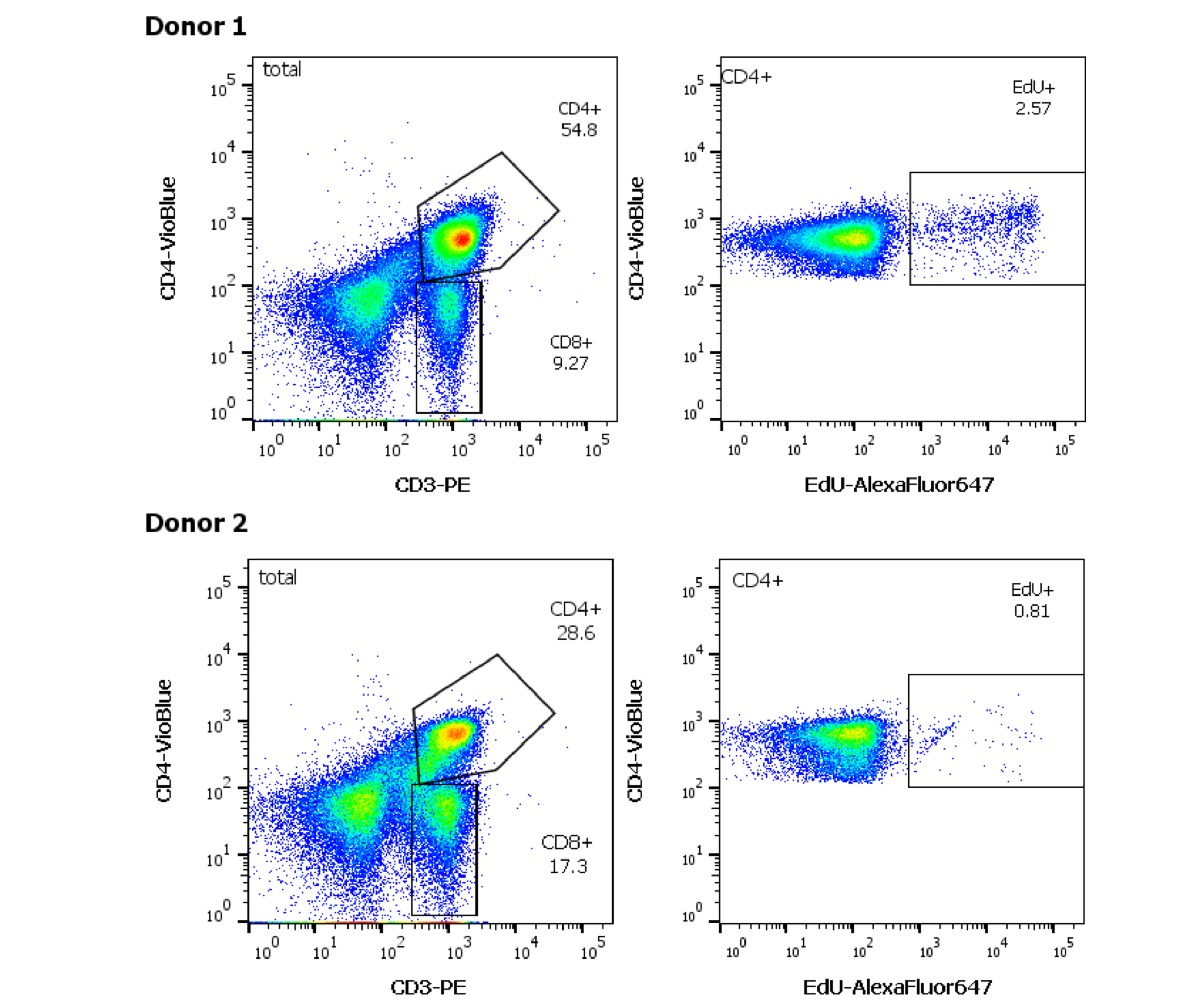
Peptide ID



- Note the highest number of peptides were identified in the AA402 – AA 421 region of the sequence.

T-Cell Activation

- Peptides presented by MHC class II are capable of stimulating T cell proliferation. One of the peptides discovered by LC-MS/MS (red box, “Peptide ID”) was synthesized and fed to PBMCs from the same donor from the discovery experiment (Donor 1) and another donor not known to present Remicade peptides (Donor 2). After 5 days in culture, T cell proliferation in the presence of this peptide was assessed by EdU incorporation analyzed by flow cytometry.



Summary

- Presented optimized methods for the identification of neoantigens (MHC class I) and biotherapeutic neopeptides (MHC class II).
- These assays have utility to improve prediction of neoantigen identification and mitigate risk of biotherapeutic immunogenicity.