

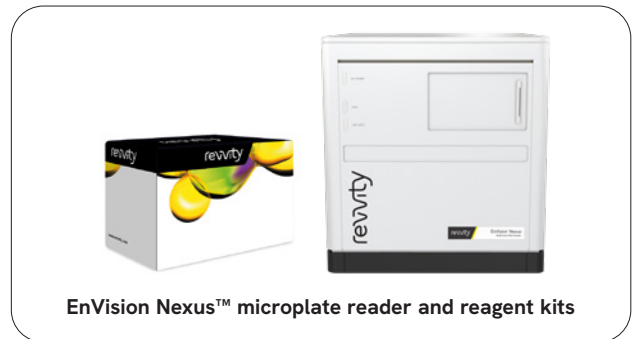


# Transforming insights into immune function.

## Solutions for immunology research

Identification of the vast array of cell types and analysis of cellular signaling molecules in the immune system is critical to the development of effective treatments for autoimmunity, allergy, and cancer. Learn about our solutions to study the pathways and network of cells.

- **BioLegend reagents:** Biomarker detection multicolor panels, enzyme-, fluorophore- and oligo-conjugates
- **BioLegend flow cytometry solutions:** Antibodies (primary and secondary) and fluorophores for panel development and phenotyping
- **Cellaca™ PLX image cytometer and reagents:** Reliable total cell and subpopulation analysis including sensitive cell counts, viability readouts, and multiplex analysis in seconds
- **Cytokine assay solutions:** Multimode microplate readers to detect signal from Alpha™, HTRF™, LANCE™, and DELFIA™ technologies including cytokine detection reagent kits
- **Fontus™ automation:** Liquid handler workstation - easy-to-use software, optimized deck access, verified protocols and reformatting capabilities to simplify your workflows and improve turnaround times



EnVision Nexus™ microplate reader and reagent kits



Cellaca PLX image cytometer



Fontus liquid handling workstation