



Accelerate T cell immunophenotyping with the Fontus and Cellaca PLX integration.

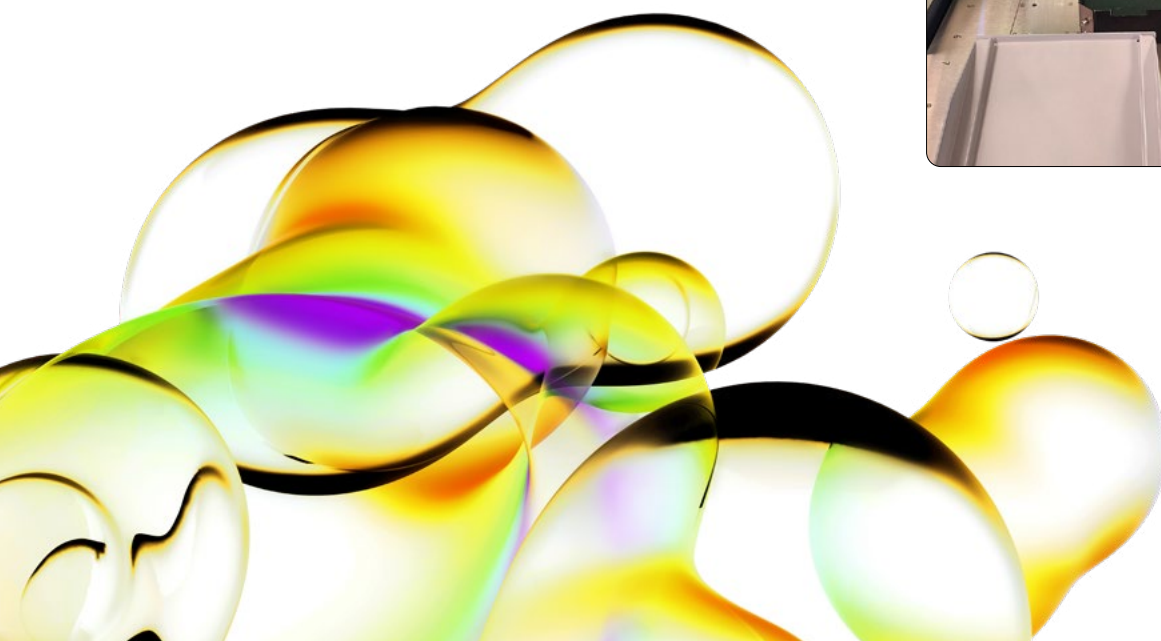
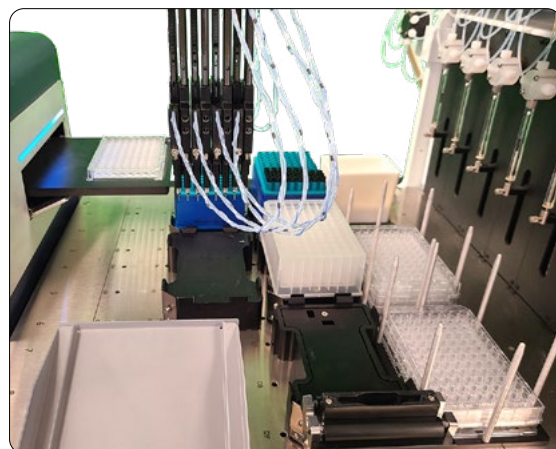
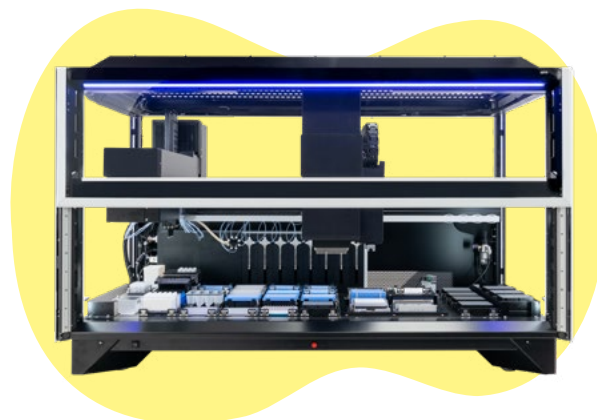
Seamless sample preparation and high-throughput analysis for cellular therapy workflows.

Transform your T cell workflow

Integrating Revvity's Fontus™ liquid handler with the Cellaca™ PLX image cytometry system enhances consistency, speed, and reproducibility. Together, they streamline immunophenotyping and viability analysis for peripheral blood mononuclear cells (PBMCs) in cell therapy research and manufacturing.

Key benefits:

- Automated staining and serial dilution of PBMCs
- Minimized hands-on time and user variability
- Fast, high-throughput imaging-based cytometry
- Comparable results to conventional flow cytometry



Automated, high-quality imaging.

Workflow overview

1. Sample prep with Fontus workstation

- Automates reagent addition, mixing, and transfers for red blood cell lysis, CD enrichment, and fluorescent antibody staining
- Loads sample to Cellaca PLX consumable for cell population analysis
- Ensures consistent timing and volume across all wells

2. Imaging and analysis with Cellaca PLX image cytometry system

- Image-based quantification and viability of CD markers for T cell, B cell, NK, Monocyte, and T cell activation markers
- Rapid processing: ~60-180 samples/hour
- Easy data export to FCS Express™ for streamlined gating and reporting

Performance

Accuracy within ~5% of leading flow cytometry platforms.
Optimized for PBMCs pre/post-RBC lysis and CD3 enrichment.

Use case: High-throughput T cell immunophenotyping

A peer-reviewed study demonstrated the use of Fontus liquid handler and the Cellaca PLX image cytometry system to perform:

- Red blood cell lysis
- CD3+ T cell enrichment (positive/negative magnetic selection)
- Multiparametric immunophenotyping (CD3, CD4, CD8)
- Viability analysis

Outcomes:

- Workflow simplified with fewer manual steps
- Throughput increased while preserving assay integrity



Why choose this integration?

Efficiency. Simplicity. Confidence.

- **Faster decisions:** move samples to downstream assays sooner
- **Lower risk:** reduce error-prone manual handling
- **Flexible scalability:** from R&D to GMP-like automation readiness

Ready to streamline your cell therapy assays?

**For more information,
contact your local
representative or
visit our website:**

revvity