

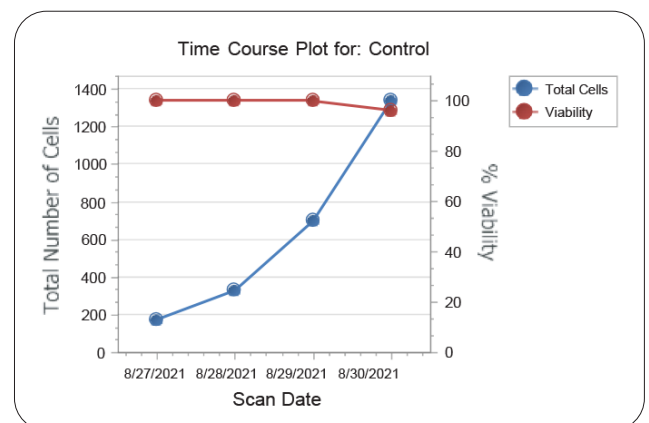


Time course graphing for improved workflow.

For the Cellaca™ MX high-throughput cell counter

Modernized software and time course graphing for improved workflow

- Effortlessly create growth curves and automatically calculate doubling time
- Measure cell concentration and viability with a few simple clicks
- Unified software platform for single-sample and high-throughput instrumentation to simplify training and documentation requirements
- Turnkey workflow solutions with scalable 21 CFR part 11 options
- Customize different counting parameters and dilution factors for every sample within a single run
- Summarize results and view custom reports to focus on actionable data for downstream work
- Streamline both low and high-throughput cell-based assays with high intra-instrument precision



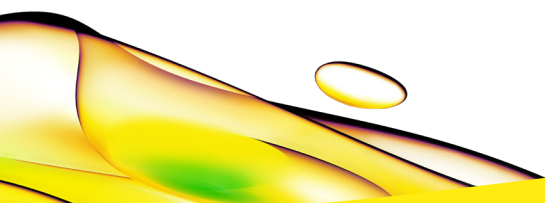


Data summary table

Growth curve tab

Elapsed time and doubling time are reported

Quick visual assessment of viability



revvity