



# The high-throughput automated cell counter that changes everything.

Count 24 samples in less than 3 minutes.

## Cellaca™ MX High-throughput automated cell counter

Your workflow will be streamlined and more efficient when you integrate the Cellaca MX with these capabilities:

- **Batch samples** - run up to 24 samples at a time
  - **Small sample volume** - only 25  $\mu$ L of cell sample required
  - **Autofocus** - fast autofocus prior to analysis
  - **Cell based assays** - data output at a single cell level enables you to do cell-based assays such as apoptosis, protein expression (including GFP and RFP), and reactive oxygen species (ROS)
  - **Multiple configurations** - choose to add up to five fluorescent filters
  - **Automation ready** - robotic integration ability with optional API
  - **Customizable reports** - add graphs, images, charts, and tables
- **Multi-language support** - over 7,000 languages available
  - **21 CFR Part 11 ready** - optional add-on that includes audit trail, user access control, and digital signature

