

CytoFLEX SRT

Fast. Easy. Accurate.

LEGACY SORTERS

CytoFLEX SRT

Power On & Instrument Warmup



20-60 minute
laser warmup

10 minute
laser warmup



QC/Sort Setup



Daily alignment of stream
or disposable cuvette*

Multiple reagents needed



Highly trained operator
verification required

20-60+ minutes

Single reagent
Fixed Optical Alignment
Single click workflow



Automated for
ease-of-use

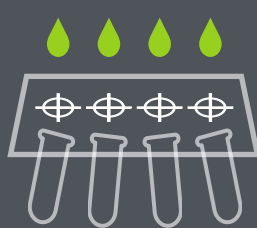


Less than 10 minutes

Ensure Stream Accuracy



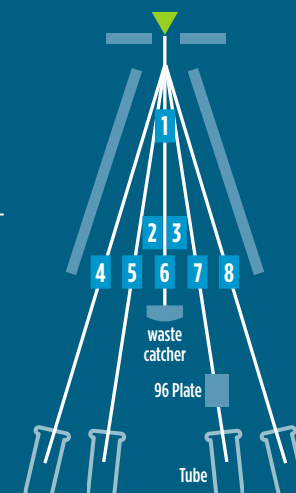
Iterative; test and fine
tune (5-20 minutes)



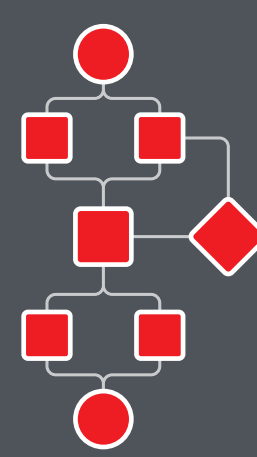
Pre-sort monitoring based
on droplet position. No
sort stream monitoring.

Automated placement
of side-streams

Side-streams and droplet
actively monitored
throughout experiment



Choosing Sort Output



Complicated
decision trees

Unintuitive,
multi-step process

Up to 384 well
plate sorting

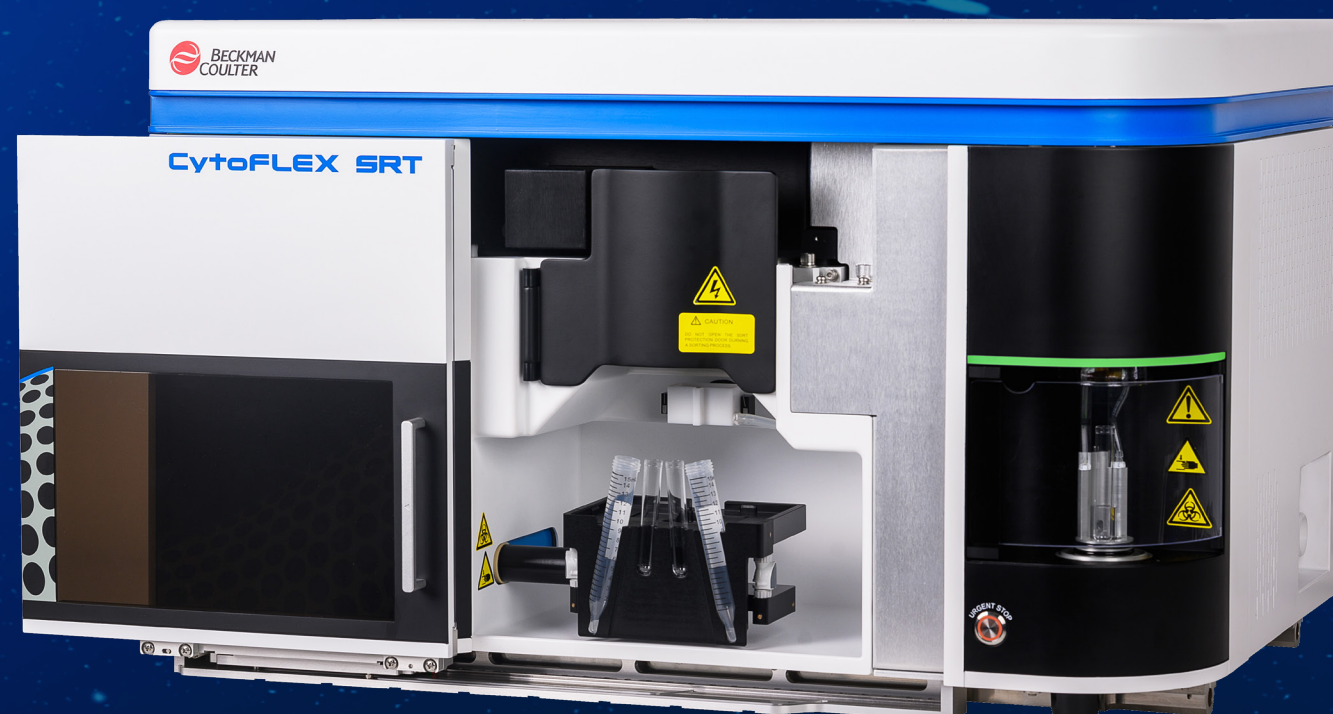
Straight down sorting
for precise deposition

Icon-based interface



Innovative Setup, Monitoring, and Stream Maintenance

CytoFLEX SRT features experience-enhancing automation that simplifies setup and operation, empowering both new and experienced researchers to focus on the experiment design without the burden of mastering a complicated instrument.



Learn more:
becl.co/cytoflex-srt

**BECKMAN
COULTER**
Life Sciences

For Research Use Only. Not for use in diagnostic procedures. *Not all legacy systems require. 21.07.2060.FLOW