

Introducing the iQue3

Get to your destination faster.



The iQue3 system is the most advanced flow cytometry platform available today with a focus on achieving actionable results in record time with an ease of use that is unparalleled. A complete platform solution of instrument, software, and reagents streamlined to rapidly generate results not just data.

Whether your destination is identifying a new drug candidate, screening for specific antibody clones, characterizing T Cells, or developing the latest immunotherapy, the iQue3 can get the answers to your research questions faster. Enabling your research to stay ahead of the curve.



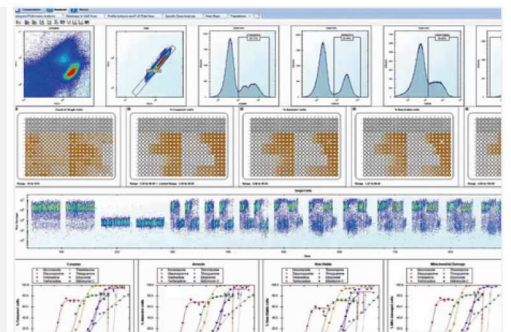
The iQue3:

- The high-throughput flow cytometry platform that combines speed and reproducibility to become the workhorse of your lab.



New! Enhanced Rinse Station:

- Up to 48 hours continuous run time.
- Reagent levels reported through the software.
- Automatic vortexing of QC beads just prior to sampling for greater consistency.
- Custom reagent cartridges reduce evaporation.



ForeCyt software (included):

- Takes data analysis from days to minutes with its powerful processing algorithms and visualization tools.

Faster, Smarter, Flow Cytometry

www.sartorius.com/intellicyt