

SARTORIUS

iQUE® 5



iQue® 5 High-Throughput Screening (HTS) by Cytometry Fluorochrome Chart

Simplifying Progress

SARTORIUS

Revealing Insight with Multiplexed Panels

Selecting the most appropriate fluorochromes for multiplexing markers of cell phenotype and function in a single well is critical to clearly resolve complex cellular research questions to support the efficient characterization of disease state.

Designing Multicolor Experiments is as Easy as 1, 2, 3:

1.

Know Your Instrument

- What are the number and type of lasers available?
- What detector | filter combinations are available to capture the signal?

This determines what labeling and detection reagents are possible to use on your system and how they will perform.

2.

Know Your Detection Reagents

- What laser wavelength excites them?
- What is the emission spectra?
- Does spillover exist between adjacent channels?
- How bright are they? What is their stain index?
- Beware that tandems and other dyes can degrade when exposed to light, temperature changes, or fixation.

This information will allow you to determine which fluorochrome will be best matched to which antigen or marker of interest.

3.

Know Your Biology

- Reserve the brightest fluorochromes for the lowest expressed antigens and vice versa.
- Avoid spillover of highly expressed antigens into detectors requiring high sensitivity.

Now, you have enough information to begin to answer your experimental questions by optimizing fluorochromes and dyes for your biology on your instrument.

This is made easier with the use of a panel builder such as the iQue® Cytometry Panel Builder (powered by EasyPanel) which is powered by intelligent algorithms that consider your instrument configuration, antigen expression levels, co-expression profiles, fluorochrome brightness, and spillover spread matrix to rapidly configure panels.



iQue® 5 HTS Platform

iQue® and Incucyte® reagents are indicated in bold

| Laser | Detector/ Filter | Channels | Fluorochromes | Fluorescent Proteins | Viability | Cell Signalling: Proliferation, Function, and Apoptosis | DNA and Cell Cycle | Incucyte® Reagents |
|-----------------|---------------------|----------|--|-----------------------------------|---|---|--|---|
| Violet 405mm | 445/45 | V445 | BD Horizon™ BV421 Pacific Blue Alexa Fluor™ 405 | eBFP Azurite Cerulean | iQue® Cell Membrane Integrity (V Blue) Dye Calcein Violet LIVE DEAD® Fixable Violet DAPI | iQue® Cell Proliferation and Encoding V Blue (Tag-it Violet™) Dye CellTrace® Violet Click-iT Plus EdUPacific Blue PolyAn Blue Multiplex Beads | DAPI FXCycle® Violet Vybrant® DyeCycle Violet | |
| | 525/45 | V525 | Pacific Green™ Horizon BV510 VioGreen® | eGFP | LIVE DEAD® Fixable Aqua | | | |
| | 586/20 | V586 | BD Horizon™ BV570 Pacific Orange | | LIVE DEAD® Fixable Yellow | | | |
| | 615/20 | V615 | BD Horizon™ BV605 Qdot™ 605 | | | | | |
| | 667/30 | V667 | BD Horizon™ BV650 Qdot™ 655 | | | | | |
| | 695/40 | V695 | Qdot™ 705 NovaFluor Violet 690 | | | | | |
| | 725/40 | V725 | BD Horizon™ BV711 Qdot™ 705 NovaFluor Violet 745 | | | | | |
| | 780/60 | V780 | BD Horizon™ BV785 Qdot™ 800 NovaFluor Violet 800 | | | | | |
| Blue 488mm | 525/45 | B525 | FITC Alexa Fluor® 488 BD Horizon™ BB515 Spark PLUS B550™ | eGFP Emerald mNeon Green | iQue® Cell Membrane Integrity (B Green) Dye Calcein AM LIVE DEAD® Fixable Green Zombie Green | iQue® Cell Proliferation and Encoding (B Green) Dye iQue® Human Caspase 3 7 Reagent iQue Qbeads® Devscreen QSav Green | Vybrant® DyeCycle Green Click-iT™ Plus EdU Alexa Fluor 488 | Incucyte® Nuclight Green |
| | 586/20 | B586 | PE eFluor 570 | eYFP | | iQue® Human Annexin V Kit iQue Qbeads® Plexscreen (detection) | | |
| | 615/20 | B615 | BD Horizon™ PE-CF594 PE Dazzle® 594 PE-Texas Red PE-Alexa Fluor 610 | | PI | | PI | |
| | 667/30 | B667 | PerCP-Cy5 PE-Cy5 eFluor PE-Alexa Fluor 647 | | | iQue® Cell Membrane Integrity (B Red) Dye | iQue® Cell Cycle Kit | |
| | 695/40 | B695 | PerCp-Vio700 PerCP-Cy5.5 PE-Cy5.5 | | 7AAD | | | |
| | 725/40 | B725 | PerCP-eFluor710 | | | | | |
| | 780/60 | B780 | PE-Cy7 PE-Vio770 | | | | | |

iQue® 5 HTS Platform

iQue® and Incucyte® reagents are indicated in bold

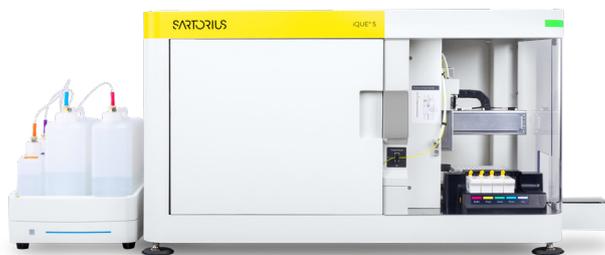
| Laser | Detector/ Filter | Channels | Fluorochromes | Fluorescent Proteins | Viability | Cell Signalling: Proliferation, Function, and Apoptosis | DNA and Cell Cycle | Incucyte® Reagents | |
|------------------|---------------------|----------|---|--|---|---|--|--|---|
| Yellow* 561nm | 586/20 | Y586 | PE Alexa Fluor® 546 | mStrawberry tdTomato AsRed 2 mRFP | CellROX® Orange | iQue® Human Annexin V Kit CellTrace Calcein Red-Orange AM | Vybrant® DyeCycle Orange Click-iT™ Plus EdU Alexa Fluor 594 | Incucyte® Nuclight Orange Incucyte® Human Fabfluor- pH Orange Dye | |
| | 615/20 | Y615 | Horizon PE-CF594 PE Dazzle 594 PE-Texas Red PE-Alexa Fluor™ 610 | mCherry HcRed mRuby mStrawberry | PI LIVE DEAD® Fixable Red | Cell Tracker Orange Cell Tracker Red | | Incucyte® Nuclight Red | |
| | 667/30 | Y667 | PE-Cy5 eFluor PE- Alexa Fluor 647 | mPlum | 7AAD | | Vybrant® DyeCycle Ruby | | |
| | 695/40 | Y695 | PE-Cy5.5 | | | | | | |
| | 725/40 | Y725 | PE-AF700 | | | | | | |
| | 780/60 | Y780 | PE-Cy7 PE-Alexa 750 PE-Vio770 | | | | | | |
| | 725/40 | V725 | Horizon BV711 Qdot 705 NovaFluor Violet 745 | | | | | | |
| | 780/60 | V780 | Horizon BV785 Qdot 800 NovaFluor Violet 800 | | | | | | |
| | Red 640nm | 667/30 | R667 | APC Alexa Fluor® 647 APC-Alexa Fluor 680 | | iQue® Cell Membrane Integrity (R Red) Dye FVD eFluor 660 LIVE DEAD® Fixable Far Red | iQue® Cell Proliferation and Encoding (R Red) Dye QuantumPlex™ StarFire Red beads (Bangs) PolyAn Red4 Multiplex beads | Click-iT™ Plus EdU Alexa Fluor 647 FxCycle® Far Red | Incucyte® Human Fabfluor- pH Red Dye |
| | | 695/40 | R695 | AF680 | | | iQue Qbeads® Plexscreen iQue Qbeads® Devscreen QSH, QSav Flash Red beads (Bangs) | DRAQ5 | |
| 725/40 | | R725 | AF700 StarBright™ Red 715 | | | | | Incucyte® Nuclight NIR | |
| 780/60 | | R780 | APC-Alexa Fluor 750 APC-Fire™ 750 APC-Cy7 APC-Vio770 StarBright™ Red 775 StarBright™ Red 815 | | FVD eFluor 780 LIVE DEAD® Fixable Near IR | iQue Qbeads® Plexscreen iQue Qbeads® Devscreen QSH | Vybrant® DyeCycle Ruby | | |

*Yellow laser is only available on 4 laser instrument version

One Platform. Unlimited Potential.

Choose the Right iQue® HTS Cytometer for Your Needs

"Rapid functional screening and immune profiling"



"A no compromise upgrade with technology and usability advances for the future of HTS using flow"

"A sweet spot in functional profiling from antibodies to CAR-T"

| | iQue® 5 (VBR) | iQue® 5 (VYBR) |
|--|--|------------------------------------|
| Ideal for complex immunology and CAR-T panel design, and for future-proofing end-to-end antibody discovery workflows | | |
| Lasers/Configuration | 3 (Violet-Blue-Red) | 4 (Violet-Yellow-Blue-Red) |
| Color Channels | 19 | 25 |
| SSC/FSC | | Yes |
| Photodetector | | SiPM |
| Particle Size Detection | | > 0.2 µm |
| Size/Dimensions (HxWxD) | | 46 x 130 x 63 cm |
| Sensitivity | | 7.2 Decades |
| Engine/Software | | Powerful Engine / Software |
| Clog Detection | | Yes, bubble sensor traces air-gaps |
| Plates | | 96-, 384-well |
| Extend Run-Time with Add-on | iQue® QMax Workstation for Extended Run-Time Workflow (Cat. Number BA-97146) | |
| Catalogue Number | BA-97134 | BA-97135 |

iQue Cytometry Panel Builder

Use the dynamic, interactive panel design tool, for streamlining the creation of complex panels.



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

Orders EMEA

Email: euorders.UK03@sartorius.com

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

Orders USA

Email: orders.US07@sartorius.com