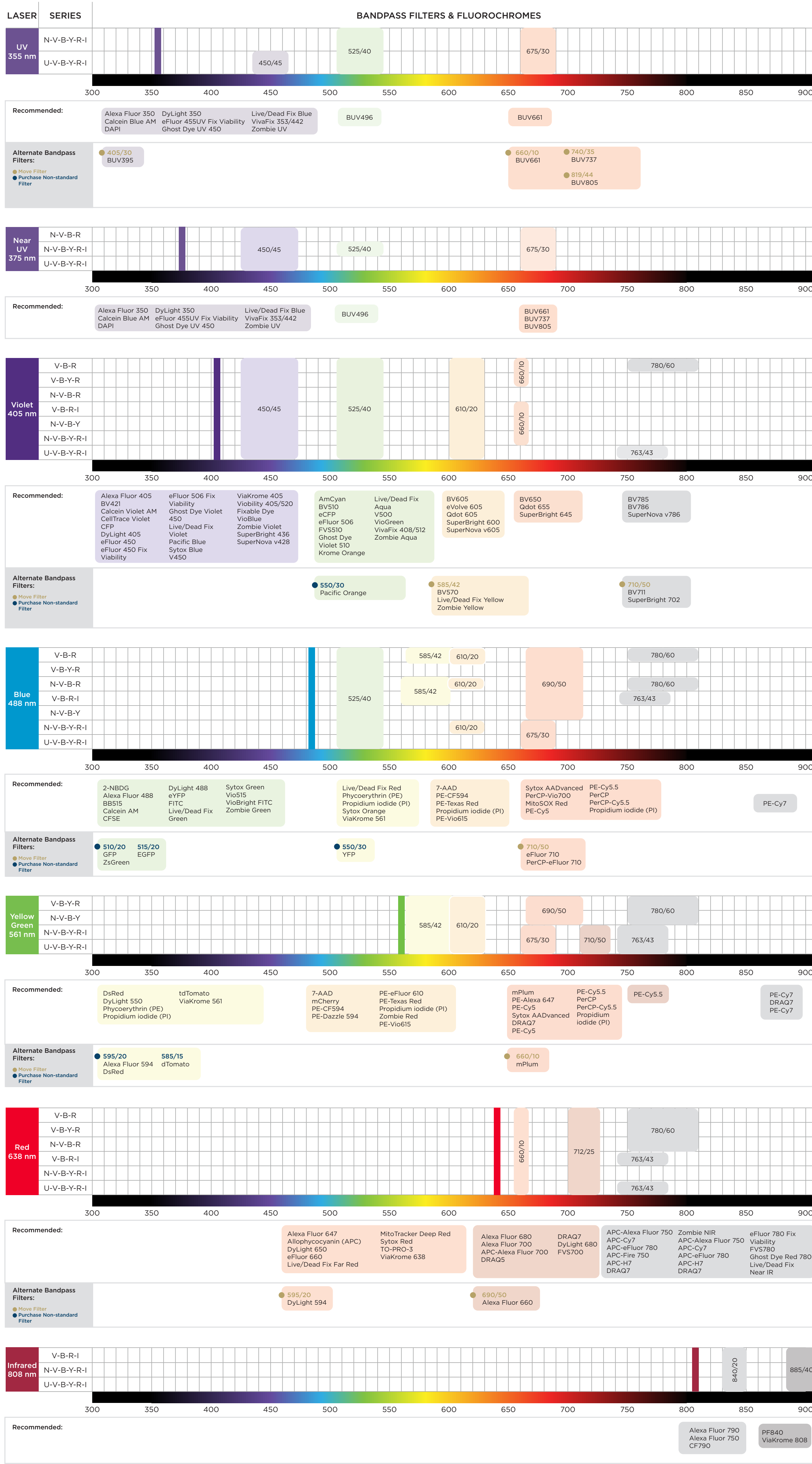


Multicolor Immunophenotyping

Panel Design Made Easy with the CytoFLEX Flow Cytometry Platform

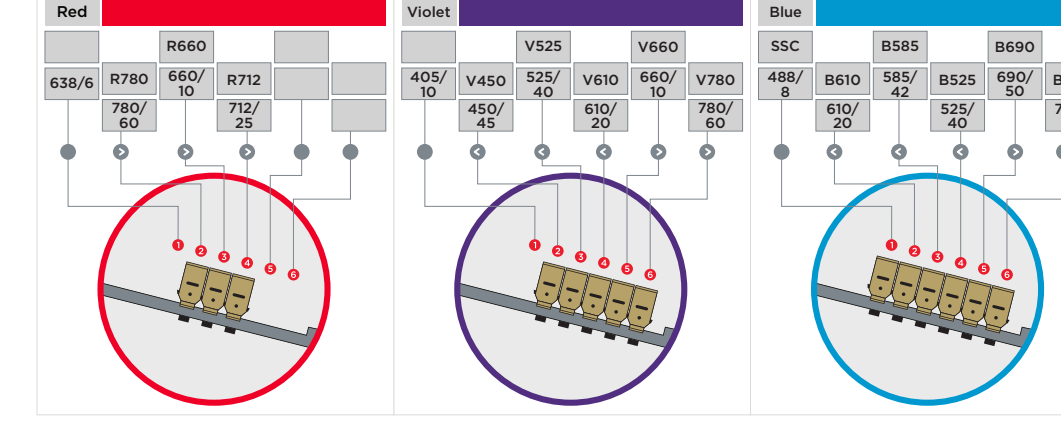
Fluorochrome Chart



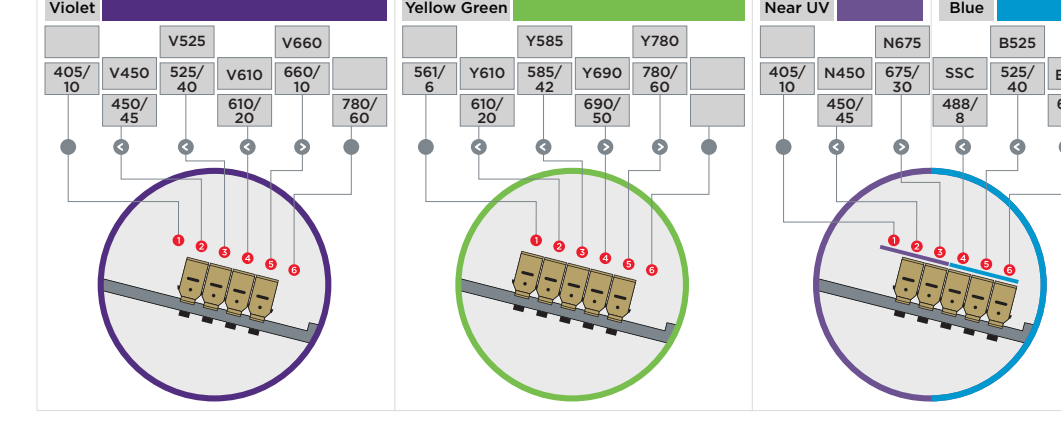
Wavelength Division Multiplex Configurations



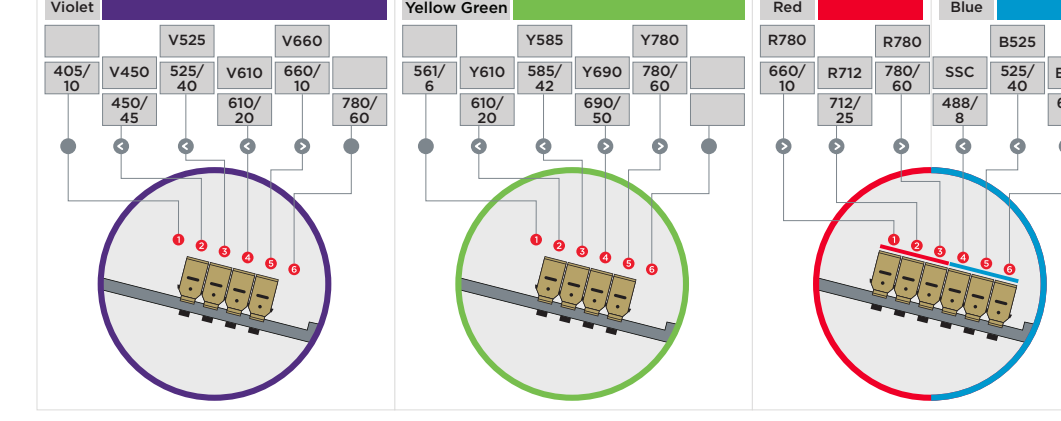
CytoFLEX Violet-Blue-Red Series



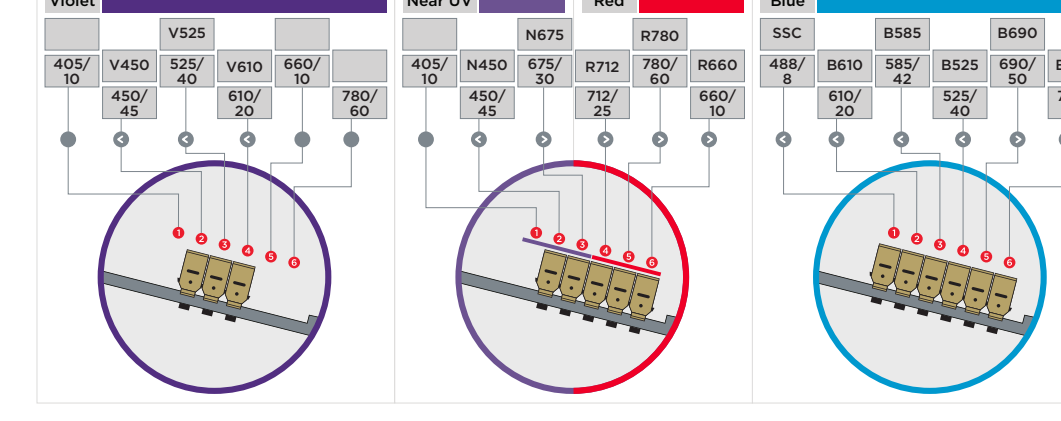
CytoFLEX S Near UV-Violet-Blue-Yellow Green Series



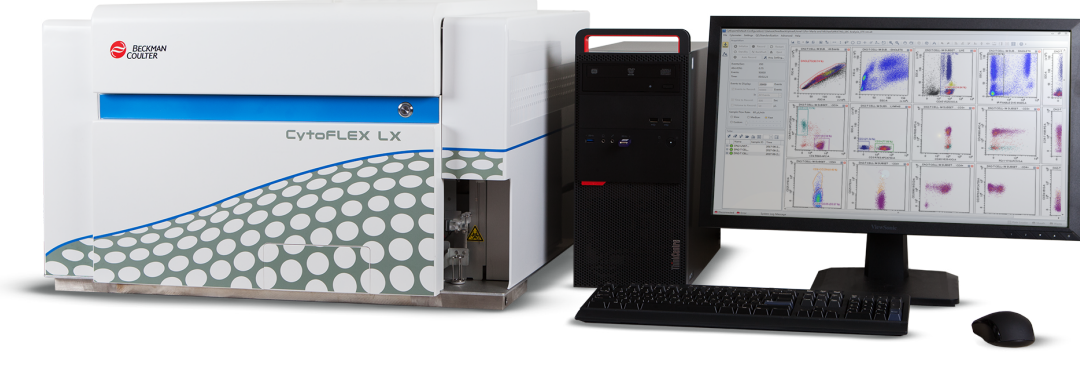
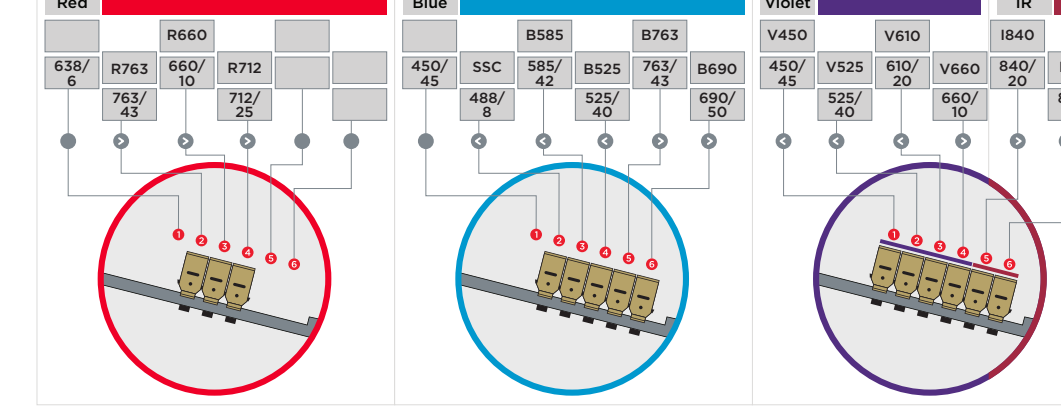
CytoFLEX S Violet-Blue-Yellow Green-Red Series



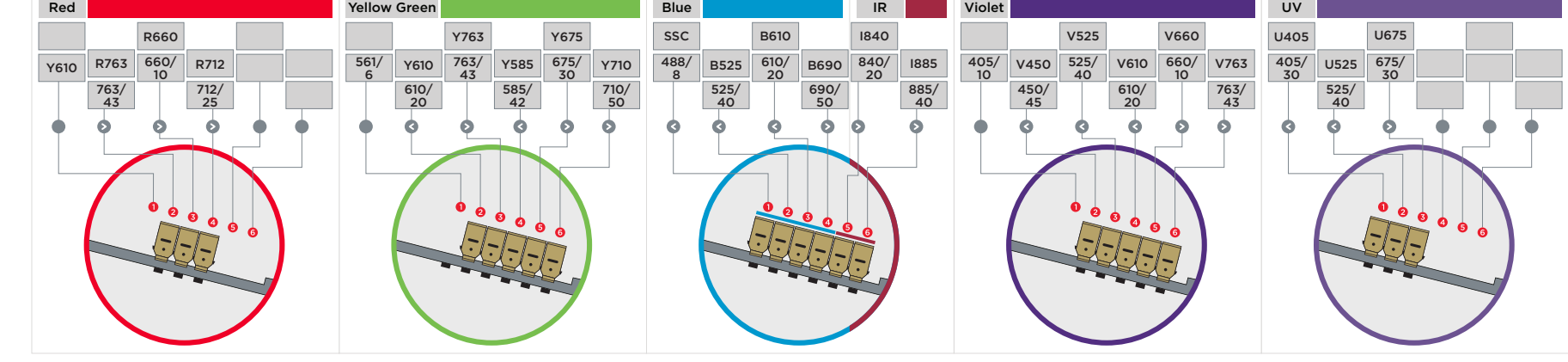
CytoFLEX S Near UV-Violet-Blue-Red Series



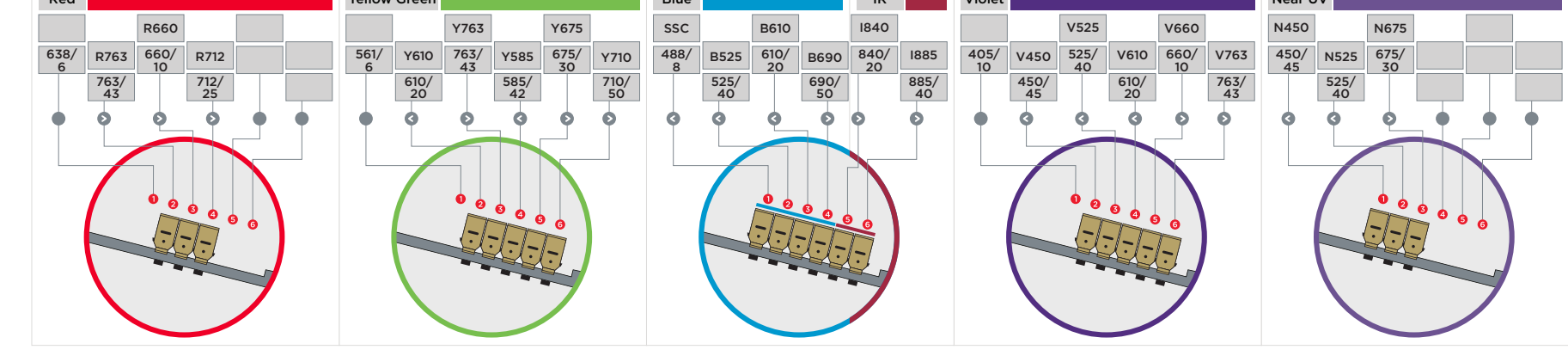
CytoFLEX S Violet-Blue-Red-Infrared Series



CytoFLEX LX UV-Violet-Blue-Yellow Green-Red-Infrared Series



CytoFLEX LX Near UV-Violet-Blue-Yellow Green-Red-Infrared Series



Bandpass Filters

Part Number	Description
A01-0048	405/10 nm Bandpass Filter
B99146	405/30 nm Bandpass Filter
A01-0050	488/8 nm Bandpass Filter
B76128	510/20 nm Bandpass Filter
B76124	515/20 nm Bandpass Filter
B76139	550/30 nm Bandpass Filter
B72627	561/6 nm Bandpass Filter
B76121	565/15 nm Bandpass Filter
B71089	585/30 nm Bandpass Filter
B76117	595/20 nm Bandpass Filter
A01-0054	638/6 nm Bandpass Filter
B78217	740/35 nm Bandpass Filter
B78220	819/44 nm Bandpass Filter
B90300	450/45 nm Bandpass Filter with Signal Attenuation (ODT)
B90294	510/20 nm Bandpass Filter with Signal Attenuation (ODT)
B90303	525/40 nm Bandpass Filter with Signal Attenuation (ODT)
B90297	610/20 nm Bandpass Filter with Signal Attenuation (ODT)
A01-0049	450/45 nm Bandpass Filter
A01-0051	525/40 nm Bandpass Filter
A01-0052	585/42 nm Bandpass Filter
A01-0053	610/20 nm Bandpass Filter
A01-0055	660/10 nm Bandpass Filter
B78244	675/30 nm Bandpass Filter
A01-0056	690/50 nm Bandpass Filter
B71092	710/50 nm Bandpass Filter
A01-0057	712/25 nm Bandpass Filter
B99143	763/43 nm Bandpass Filter
A01-0058	780/60 nm Bandpass Filter
B99144	840/20 nm Bandpass Filter
B99145	885/40 nm Bandpass Filter
C30171	Custom Optical Filter Holder (1) with Screws (2)
C30249	Custom Optical Filter Holder Mounting Fixture
C32857	Custom Optical Filter Holder Screws (2)
C32857	Custom Optical Filter Holder Screws (2)

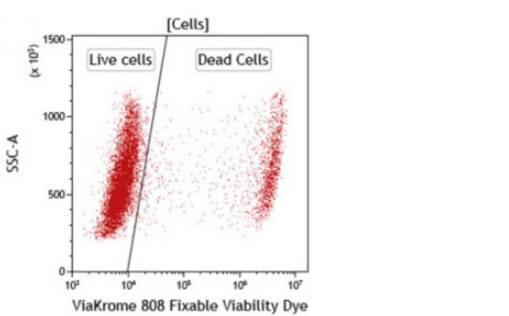
Custom Filter Holder

Part Number	Description
C30171	Custom Optical Filter Holder (1) with Screws (2)
C30249	Custom Optical Filter Holder Mounting Fixture
C32857	Custom Optical Filter Holder Screws (2)

ViaKrome Fixable Viability Dyes

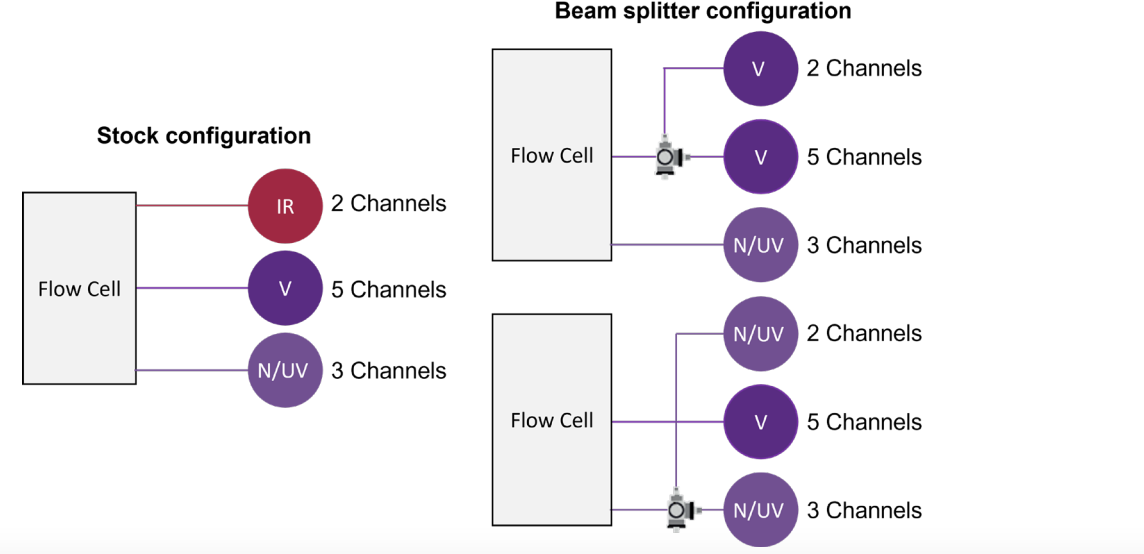
- Compatible with surface and intracellular staining
- Improved performance over standard NHS ester-based dyes
- Soluble in biologically compatible buffer, PBS
- Does not require organic solvents such as DMSO for reconstitution

beckman.com/cell-health



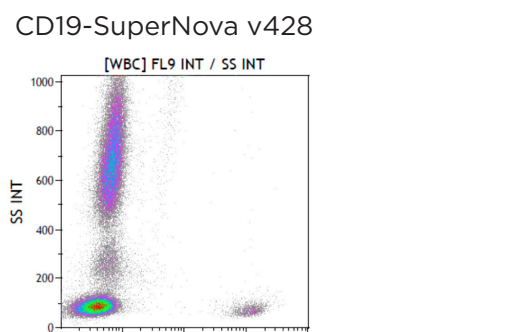
Expand Violet, UV or Near UV channels with the CytoFLEX LX Beam Splitter

- Addition of the CytoFLEX LX Beam Splitter allows the instrument to configure the IR detectors to detect emission from the UV, Near UV, or Violet WDM
- User can switch between extending the violet or the N/UV channels



SuperNova Dyes

- Deliver the highest staining index on the violet laser for their respective channel
- Provide greater confidence in results thanks to a proprietary formulation minimizing nonspecific staining
- Available as catalog products or can be provided within dry custom multicolor panels



CytoFLEX SRT Cell Sorter Violet-Blue-Yellow Green-Red Series

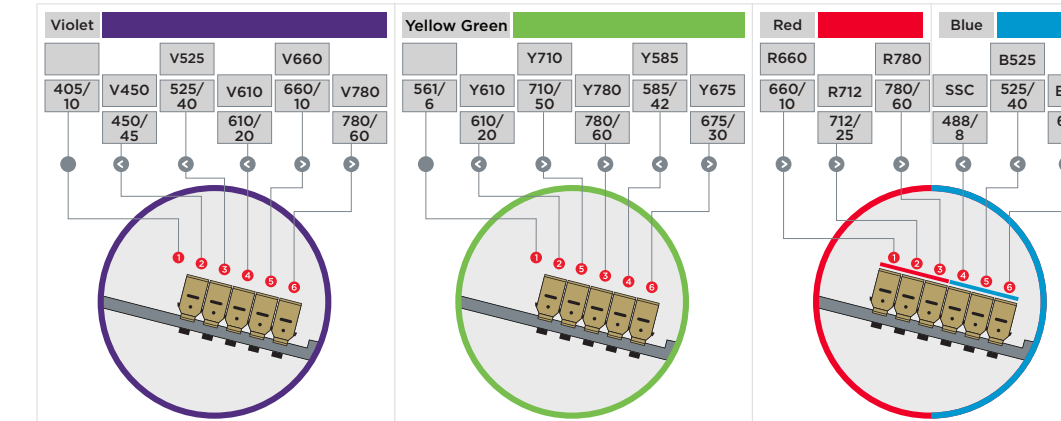


DIAGRAM KEY

- Use for bandpass filters with wavelengths shorter than 660 nm
- Use for bandpass filters with wavelengths longer than 660 nm
- Use for spare bandpass filter storage

Instrument Quick Start Guide

Startup	Verify System Performance	Create Compensation	Set up Experiment	Shut Down
<ol style="list-style-type: none"> Turn on Start the computer Check fluid levels Select Cytometer>System Startup Program 8 minutes for system priming 	<ol style="list-style-type: none"> QC/Standardization>Start Select Lot Num>Start File>Close QC/Standardization 1-2 minutes 	<ol style="list-style-type: none"> File>New Compensation Rename experiment (if desired)>Save (.xlt) Compensation Setup >select needed channels and specify sample type>OK Initialize (if standby state) Add sample, Select the appropriate Tube>Run (if in Ready state) After acquiring all tubes, Settings>Compensation Calculation Save to Compensation Library (for future experiments) or Save as (for exporting to third party software) 	<ol style="list-style-type: none"> File>New Experiment Rename experiment (if desired)>Save (.xlt) Settings>Set Channel>add Labels as needed, deselect unused channels if desired Settings>Compensation Matrix (OR comp matrix icon from tube screen)>Choose appropriate Option>OK Create plots Add tube Establish stop criteria, set sample flow rate, desired gains (if not imported with compensation) and threshold 	<ol style="list-style-type: none"> Prepare FlowClean (2 mL) and DI water tubes (2 mL) Cytometer>Daily Clean Adjust Duration as needed Add water when instructed (step 2) When finished, Close CytoFLEX LX: Cytometer>turn off CytoFLEX or CytoFLEX S: Exit software and manually shut down instrument



Please note: The channel naming convention is based on CyExpert v2.5.
 For Research Use Only. Not for use in diagnostic procedures.
 © 2023 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.
 For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
 2022-GBL-EN-100218-v2