

# CF<sup>®</sup> Dyes for Flow Cytometry

This chart serves as a helpful reference guide for building flow cytometry panels using our bright CF<sup>®</sup> dyes and other commonly used fluorophores. We have included common laser lines and detection channels for reference.

## Emission Filters

- Pacific Blue™ 450/50
- FITC, AmCyan 525/50
- PE 575/26
- Texas Red® 610/20
- Cy5, APC, PerCP 660/20
- Cy7 780/60
- IR840 840/20
- IR885 885/40

## Laser Lines

- ✦ UV: 355 nm
- ✦ Violet: 405 nm
- ✦ Blue: 488 nm
- ✦ Yellow: 561 nm
- ✦ Red: 640 nm
- ✦ Infrared: 808 nm
- ✦ Yellow/Green: 532 nm

