



Stain Index for StarBright UltraViolet Dyes

Quick Guide

StarBright Dyes are new, proprietary fluorescent nanoparticles conjugated to Bio-Rad's highly validated antibodies. Designed with exceptional brightness, narrow excitation and emission, stability, and with high lot-to-lot reproducibility.

This handy reference guide shows how the stain index of StarBright UltraViolet Dyes compares to other ultraviolet laser-excitable fluorescent dyes that have similar maximal emissions.

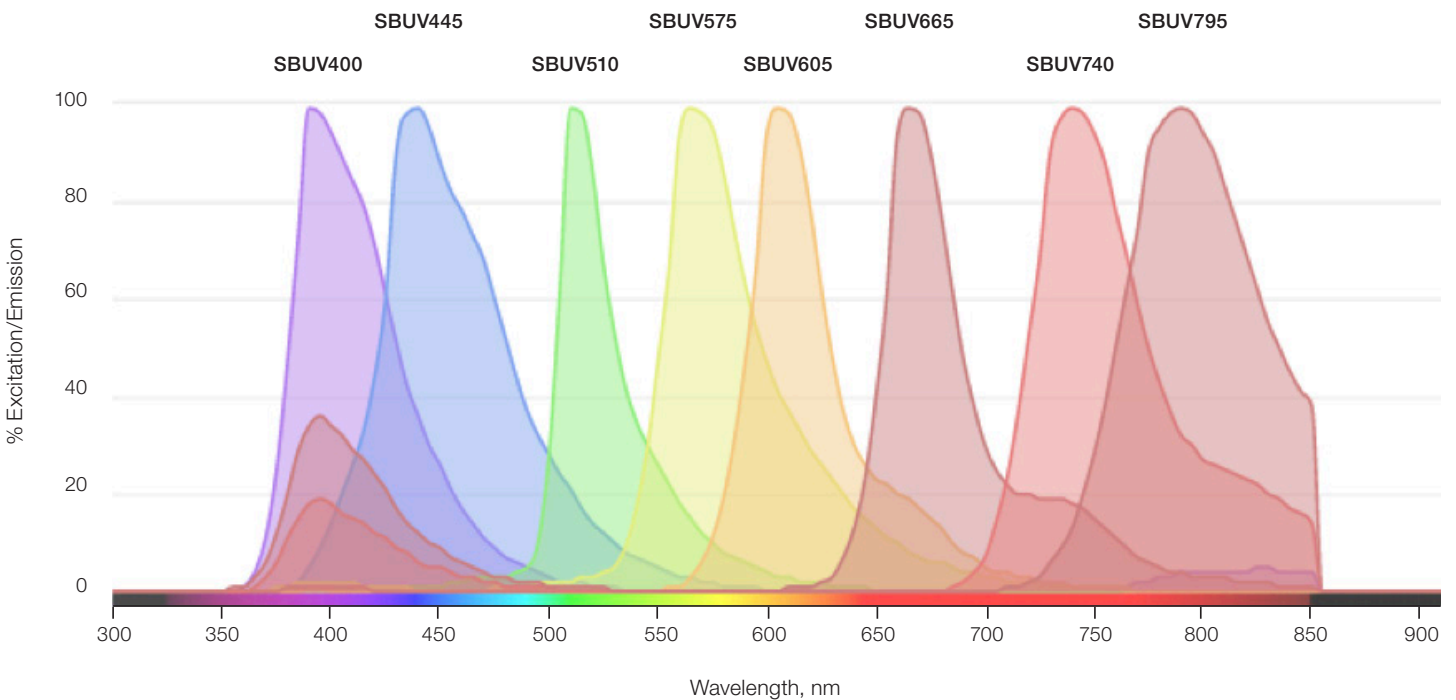


Fig. 1. Excitation and emission spectra of StarBright UltraViolet Dyes.

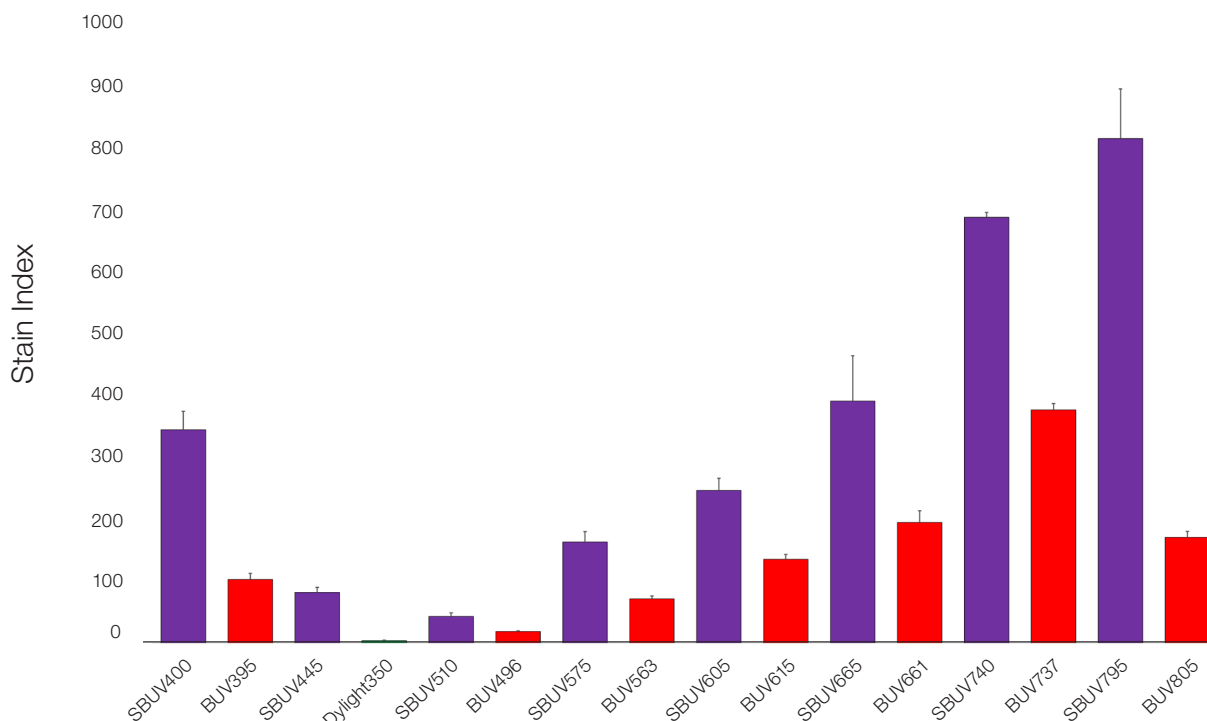
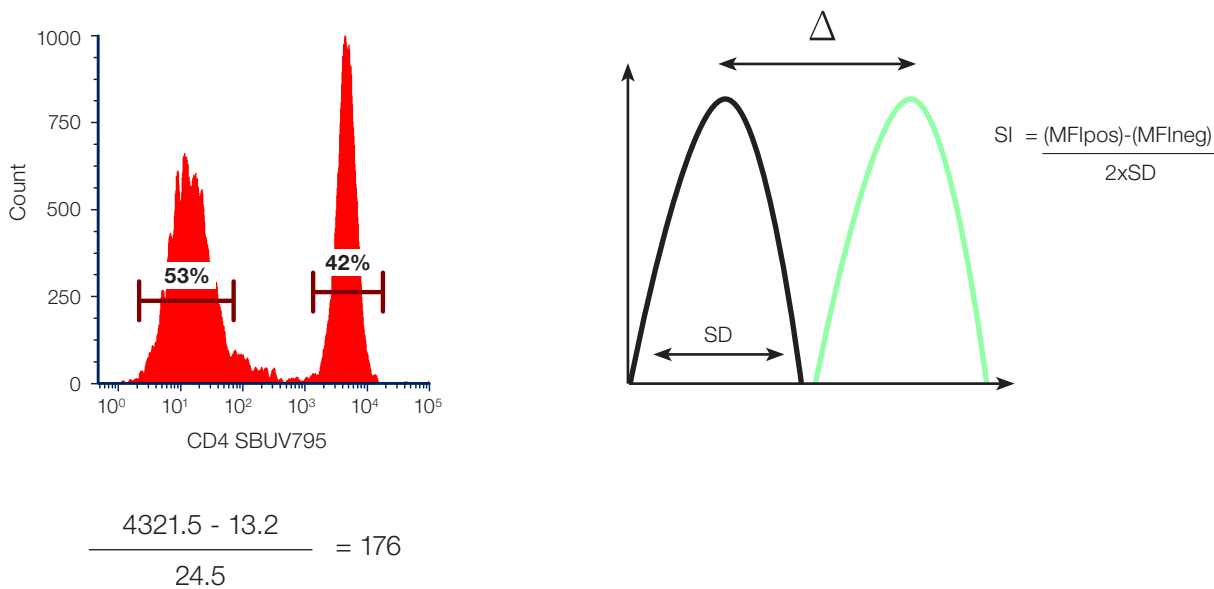


Fig. 2. Stain index comparison. Stain index of human peripheral blood stained with various 355 nm excitable StarBright UltraViolet, Brilliant Ultraviolet, and conventional dyes conjugated to CD4 (MCA1267). Data generated are the average of 3–4 donors, antibodies were titrated before use and analyzed on the ZE5 Cell Analyzer. Purple, StarBright Dyes; red, Brilliant Violet; green, other non-polymer dyes. SBUV, StarBright UltraViolet Dye; BUV, Brilliant Ultraviolet.

How to generate a stain index



Visit [bio-rad-antibodies.com/stain-index](https://www.bio-rad-antibodies.com/stain-index) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**