



Stain Index for StarBright Blue Dyes

Quick Guide



StarBright™ Dyes are new, proprietary fluorescent nanoparticles conjugated to the Bio-Rad™ line of highly validated antibodies. StarBright Dye conjugated antibodies are designed to offer exceptional brightness, narrow excitation and emission, and stability with high lot-to-lot reproducibility.

This handy reference guide shows the stain index of StarBright Blue Dyes (SBB) compared to other blue laser-excitable fluorescent dyes with similar maximal emissions.

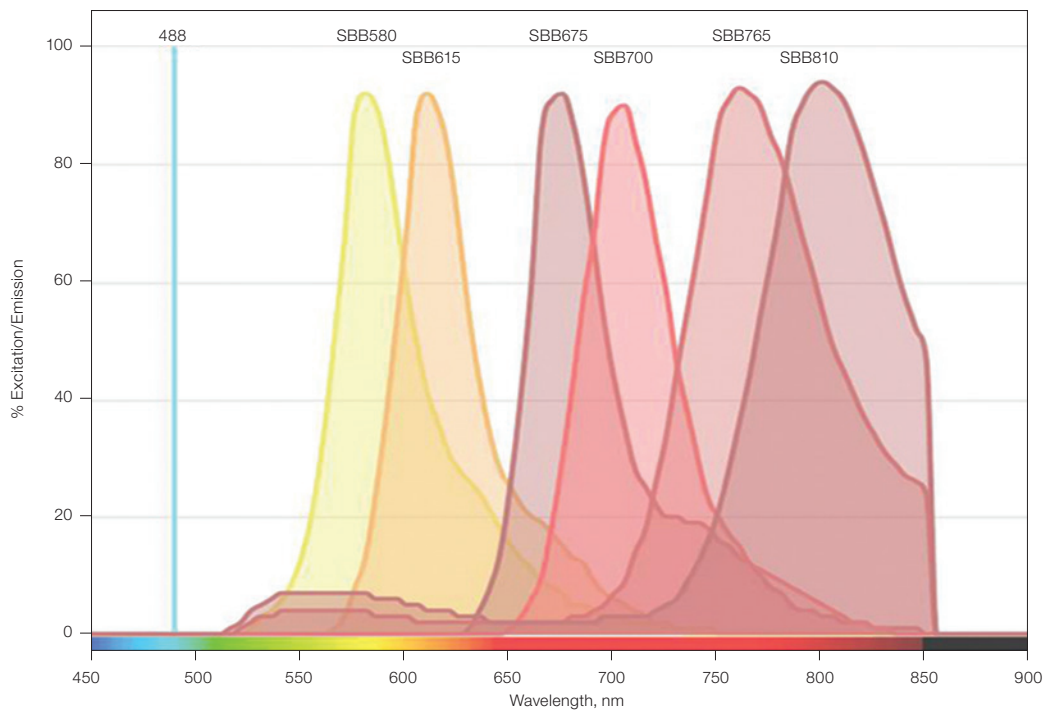


Fig. 1. Excitation and emission spectra of StarBright Blue Dyes.

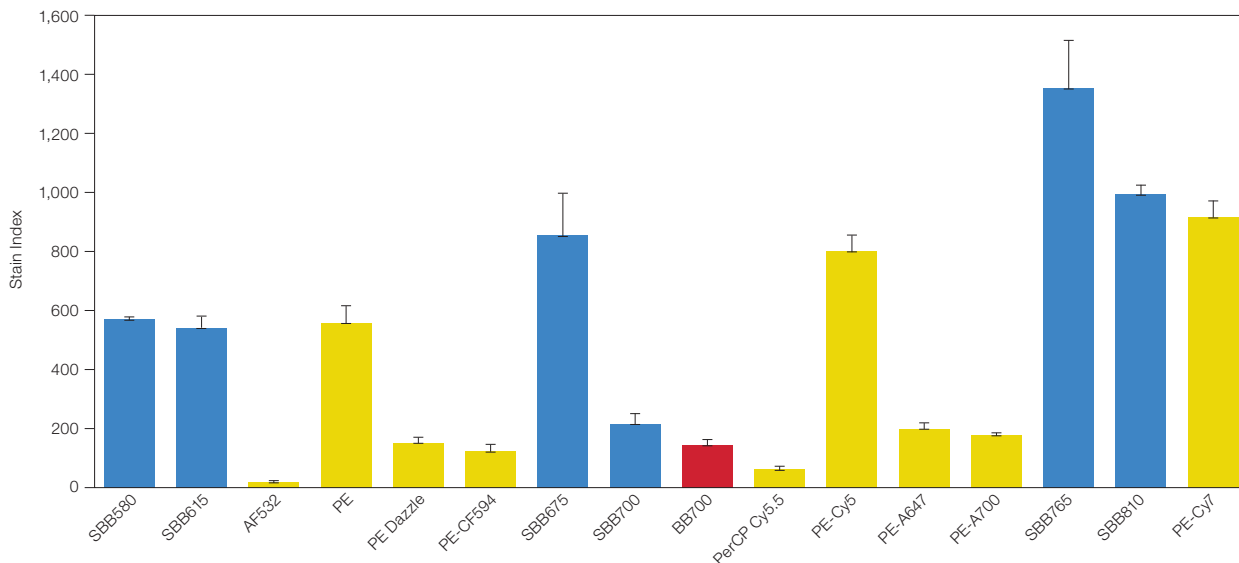
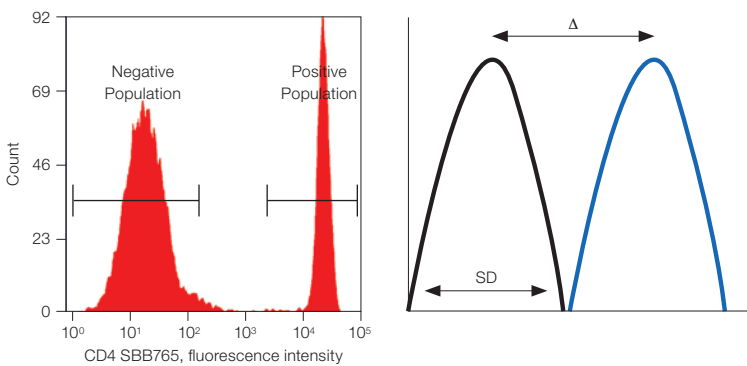


Fig. 2. Stain index comparison of various 488 nm excitable dyes used to detect CD4 in human peripheral blood samples. StarBright Blue Dyes (■), Brilliant Blue (■), and conventional nonpolymer dyes (■) were conjugated to Mouse Anti-Human CD4 Antibody (Bio-Rad Laboratories, Inc., #MCA1267). Data shown are the average of three donors. Antibodies were titrated prior to use and analyzed on the ZE5 Cell Analyzer. AF, Alexa Fluor; BB, Brilliant Blue; Cy, cyanine dye; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; SBB, StarBright Blue Dye.

How to Generate a Stain Index (SI)

A stain index is a method to calculate the relative brightness of a fluorescent dye on a particular instrument. It is defined as the difference (delta) between the mean fluorescent intensity (MFI) of the positive population (blue) and the negative population (black), divided by two times the standard deviation (SD) of the negative population. A titration of antibody followed by the SI calculation can determine the optimal amount of antibody to use in your experiments.



$$SI = \frac{(MFI_{pos}) - (MFI_{neg})}{2 \times SD_{neg}} = \frac{17,923.8 - 10.4}{2 \times 14.9} = 601$$

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

BIO-RAD and STARBRIGHT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

