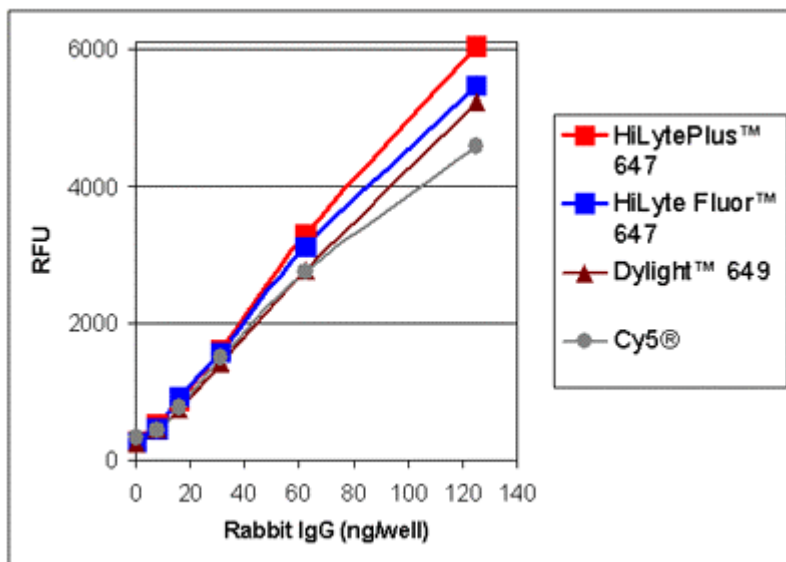




HiLytePlus(tm) Long Wavelength Dyes - Brilliant Gets Better

In head to head comparisons between HiLyte Fluor(tm) dyes and comparable dyes such as Cy[®] & others, HiLyte Fluor(tm) dyes have delivered superior results that can best be described as brilliant. With the new HiLytePlus(tm) series of long wavelength dyes, brilliant gets even better.



AnaSpec is pleased to introduce **HiLytePlus(tm)**. Built on HiLyte Fluor(tm) technology, the HiLytePlus(tm) series represents AnaSpec's highest performing dyes ever for long wavelengths.

HiLytePlus™ Series

- Even brighter than HiLyte Fluor(tm) dyes
- Like HiLyte Fluor(tm) series, outperforms Cy[®] & DyLight(tm) dyes

- Superior photostability
- Excitation wavelengths of 555, 647 and 750 nm
- Ideal for high-performance biomarker labeling
- Available as individual dyes, AnaTag(tm) Labeling Kits, labeled secondary antibodies, and labeled streptavidin.

