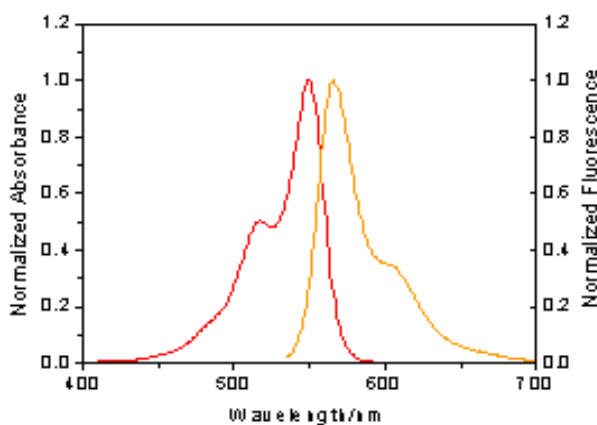




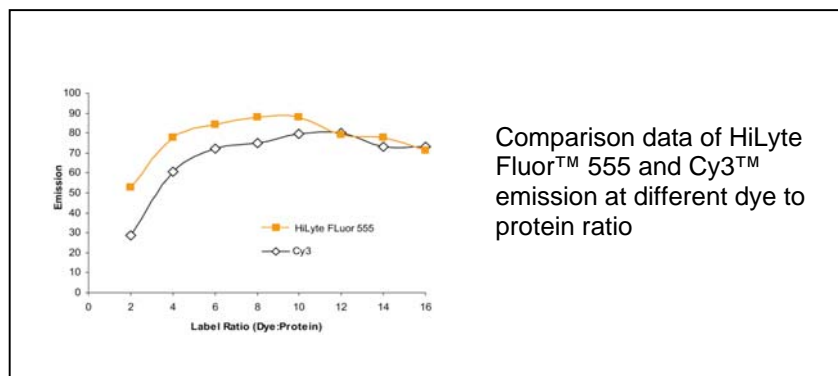
## Product Data Sheet

---

Product Name:	HiLyte Fluor™ 555 hydrazide
Size:	1 mg
Catalog Number:	81253
Molecular Weight:	883.13
Solvent:	Water or DMF
Product Description:	HiLyte Fluor™ 555 hydrazide is a carbonyl-reactive fluorescent labeling dye that can be used to generate conjugates that are slightly red-shifted compared to those of Cy3™ dyes, resulting in optimal match to filters designed for Cy3™ dyes. It can be used for labeling glycoproteins such as HRP. HiLyte Fluor™ 555 dye is more photostable than Cy3™, providing researchers with additional time for image capture. Additionally, in-house data indicated that HiLyte Fluor™ 555 is superior to Cy3™ in fluorescence polarization-based assays.
Spectra:	Maximum Ex/Em wavelength = 551/567 nm



**Flourescence Intensity:** Performance of dye-goat anti-rabbit IgG conjugates



**Stock Solution Preparation:** Prepare 10 mg/ml dye stock solution in water or DMF. Completely dissolve all dye content by vortexing. Dye solution should be prepared fresh, immediately before an experiment. Long-term storage in solution form may reduce dye activity. Any solution containing the dye should be protected from light.

**Shelf Life and Storage:** Shelf life is two years if stored at -20C, desiccated and protected from light.

**Related Products:**

Catalog#	Product Name
81250	HiLyte Fluor™ 555 acid
81251	HiLyte Fluor™ 555 acid, SE
81252	HiLyte Fluor™ 555 amine
81254	HiLyte Fluor™ 555 C2 maleimide

This product is for *in vitro* research use only