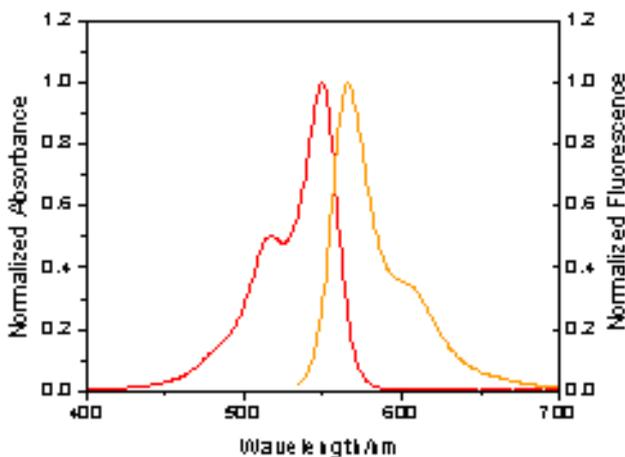




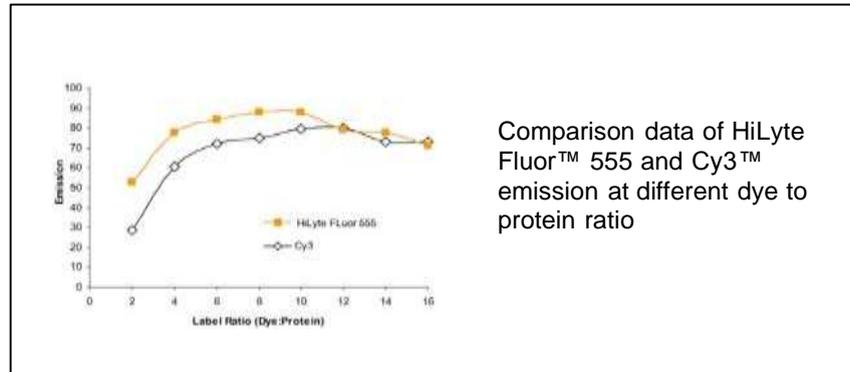
## Product Data Sheet

---

Product Name:	HiLyte Fluor™ 555 C2 maleimide
Size:	1 mg
Catalog Number:	AS-81254
Molecular Weight:	1062.3
Solvent:	Water or DMF
Product Description:	HiLyte Fluor™ 555 C2 maleimide is an excellent thiol-reactive fluorescent labeling dye that can be used to generate protein conjugates that are slightly red-shifted compared to those of Cy3™ dyes, resulting in optimal match to filters designed for Cy3™ dyes. HiLyte Fluor™ 555 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 = 552/569 nm



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



Stock Solution Preparation:

Prepare 10 mg/ml dye stock solution in high quality DMF or water. 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
AS-81250	HiLyte Fluor™ 555 acid
AS-81251	HiLyte Fluor™ 555 acid, SE
AS-81252	HiLyte Fluor™ 555 amine

This product is for *in vitro* research use only.