

## **Product Data Sheet**

Product Name: HiLyte Fluor™ 555 acid

Size: 5 mg

Catalog Number: AS-81250

Molecular Weight: 869.1

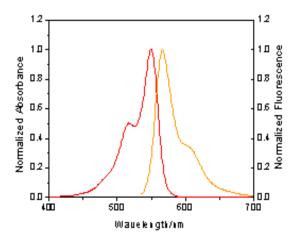
Solvent: Water

Product Description: Spectra of HiLyte Fluor™ 555 conjugates are slightly red-shifted

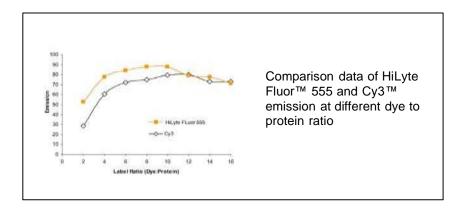
compared to those of Cy3<sup>™</sup> dyes, resulting in optimal match to filters designed for Cy3<sup>™</sup> dyes. HiLyte Fluor<sup>™</sup> 555 is more photostable than Cy3<sup>™</sup>, providing researchers with additional time for image capture. Additionally, in-house data indicated that HiLyte Fluor<sup>™</sup> 555 is superior

to Cy3™ in fluorescence polarization-based assays.

Spectra: Maximum Ex/Em wavelength = 550/566nm



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



Stock Solution Preparation: Prepare 10 mg/ml dye stock solution in 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-81251	HiLyte Fluor™ 555 acid, SE
AS-81252	HiLyte Fluor™ 555 amine
AS-81254	HiLyte Fluor™ 555 C2 maleimide

This product is for *in vitro* research use only