

Product Data Sheet

Product Name: HiLyte Fluor™ 555 amine

Size: 1 mg

Catalog Number: AS-81252

Molecular Weight: 911.18

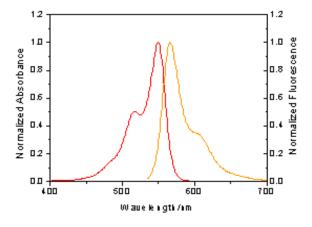
Solvent: Water

Product Description: HiLyte Fluor™ 555 amine is a carbonyl-reactive fluorescent labeling dye

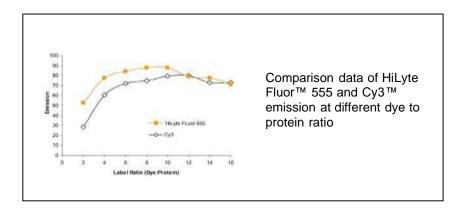
that can be used to generate conjugates that are slightly red-shifted compared to those of Cy3 dye, resulting in optimal match to filters designed for Cy3 dyes. HiLyte Fluor™ 555 is more photostable than Cy3™, providing researchers with additional time for image capture. Additionally, in-house data indicated that HiLyte Fluor™ 555 is superior

to Cy3™ in fluorescence polarization-based assays.

Spectra: Maximum Ex/Em wavelength = 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 high quality anhydrous DMSO. 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-81254	HiLyte Fluor™ 555 C2 maleimide

This product is for in vitro research use only