



Product Data Sheet

Product Name: HiLyte Fluor™ 488 C2 maleimide

Size: 1 mg

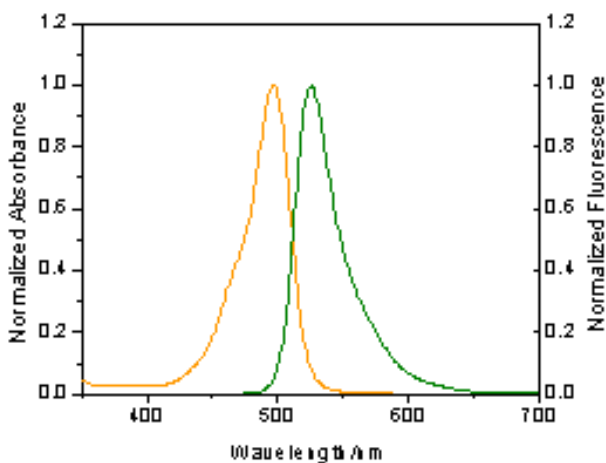
Catalog Number: AS-81164

Molecular Weight: 567.55

Solvent: DMF or DMSO

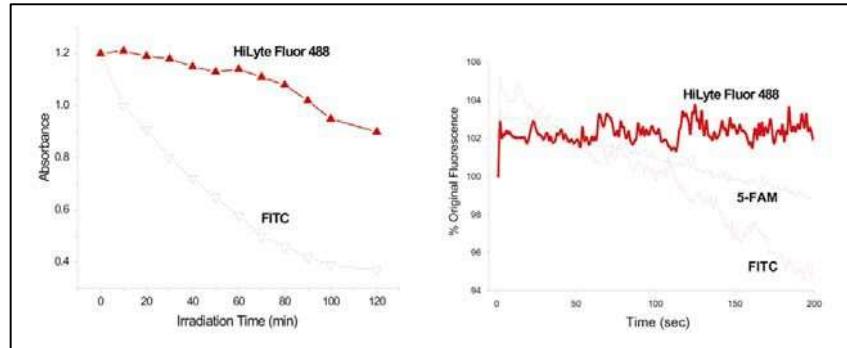
Product Description: HiLyte Fluor™ 488 C2 maleimide is an excellent thiol-reactive fluorescent labeling dye that generates protein conjugates far superior than those of fluorescein derivatives such as FITC. Not only are HiLyte Fluor™ 488 conjugates significantly brighter than fluorescein conjugates, they are much more photostable. Also, fluorescence of HiLyte Fluor™ 488 fluorophore is not pH dependent at pH range between 4 to 10. This pH insensitivity is a major improvement over fluorescein, which emits its maximum fluorescence only above pH 10.

Spectra: Maximum Ex/Em wavelength = 502/527 nm



Photostability:

HiLyte Fluor™ 488, 5-FAM and FITC photostability over time



Stock Solution Preparation:

Prepare 10 mg/ml dye stock solution in 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-81160	HiLyte Fluor™ 488 acid
AS-81161-1 (1 mg)	HiLyte Fluor™ 488 acid,SE
AS-81161 (5 mg)	HiLyte Fluor™ 488 acid,SE
AS-81162	HiLyte Fluor™ 488 amine, TFA salt
AS-81163	HiLyte Fluor™ 488 hydrazide
AS-81164	HiLyte Fluor™ 488 C2 maleimide

This product is for *in vitro* research use only