

Product Data Sheet

- Product Name: HiLyte Fluor™ 488 amine, TFA Salt
- Size: 1 mg

Catalog Number: AS-81162

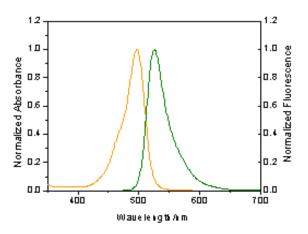
Molecular Weight: 530.45

Solvent: DMF or DMSO

Product Description: HiLyte Fluor[™] 488 amine is a carbonyl-reactive fluorescent labeling dye. Fluorescence of HiLyte Fluor[™] 488 fluorophore is not pH dependent at pH range between 4-10. This pH insensitivity is a major improvement over fluorescein, which emits its maximum fluorescence only above pH 9.

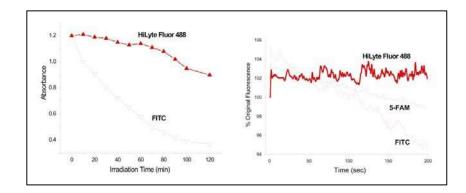
Spectra:

Maximum Ex/Em wavelength = 503/528 nm



Photostability:

HiLyte $\mathsf{Fluor}^{^{\mathrm{\tiny{TM}}}}$ 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.

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

Related Products:

This product is for in vitro research use only