

Product Data Sheet

Product Name: HiLyte Fluor™ 488 hydrazide

Size: 1 mg

Catalog Number: AS-81163

Molecular Weight: 388.38

Solvent: DMF or DMSO

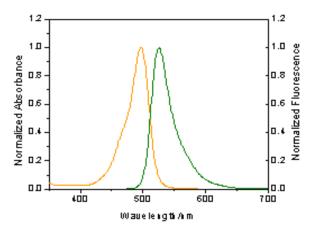
Product Description: HiLyte Fluor™ 488 hydrazide is a carbonyl-reactive fluorescent labeling

dye. It can be used for labeling glycoproteins such as HRP.

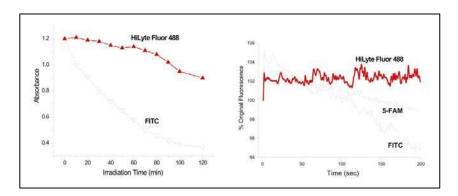
Fluorescence of HiLyte Fluor[™] 488 fluorophore is not pH dependent at pH range between 4 to10. This pH insensitivity is a major improvement over fluorescein, which emits its maximum fluorescence only above pH

9.

Spectra: Maximum Ex/Em wavelength is 503/528 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