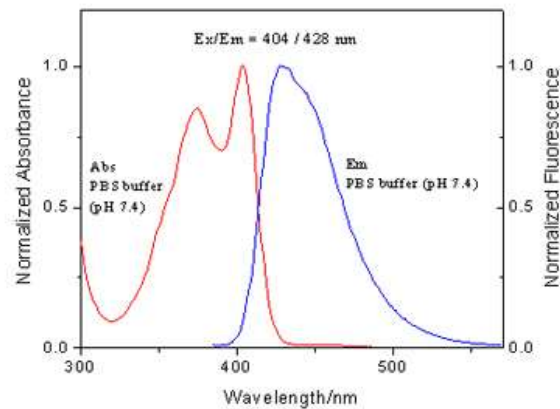




## Product Data Sheet

---

Product Name:	HiLyte Fluor™ 405 hydrazide
Size:	1 mg
Catalog Number:	AS-89319
Molecular Weight:	636.63
Solvent:	DMF or DMSO
Product Description:	HiLyte Fluor™ 405 is a blue fluorescent dye with Ex/Em = 404/428 nm, spectrally similar to Alexa Fluor™ 405 and DyLight Fluor™ 405 dyes. Conjugates of HiLyte Fluor™ 405 are well excited by the 407 nm spectral line of the Krypton laser and the 408 nm line of the violet laser diode. HiLyte Fluor™ 405 hydrazide is a carbonyl-reactive fluorescent labeling dye. It can be used for labeling glycoproteins such as HRP.
Spectra:	Maximum Ex/Em wavelength is 404/428 nm.



**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-89317-1 (1 mg)	HiLyte Fluor™ 405 acid, SE

This product is for *in vitro* research use only