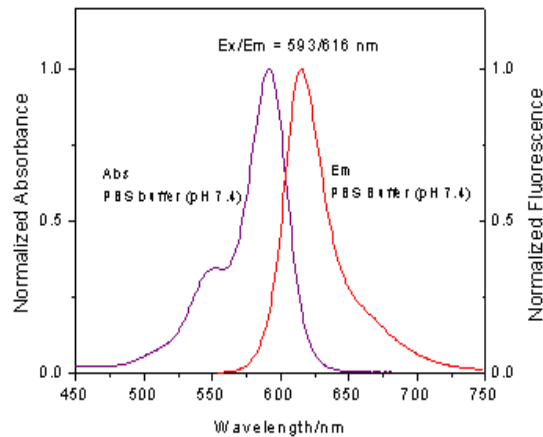




## Product Data Sheet

---

Product Name:	HiLyte Fluor™ 594 hydrazide
Size:	1 mg
Catalog Number:	AS-81274
Molecular Weight:	879.9
Solvent:	DMF or DMSO
Product Description:	HiLyte Fluor™ 594, an alternative to Alexa Fluor™ 594 and DyLight Fluor™ 594, has spectral characteristics similar to those of Texas Red. Its labeling performance and stability are better than those of Texas Red. It has high extinction coefficient (80,000 M <sup>-1</sup> cm <sup>-1</sup> ), high fluorescence quantum yield (0.9) and low correction factor (0.17). HiLyte Fluor™ 594 based bioconjugates exhibit little spectral overlap with green fluorescent conjugates, and can be efficiently excited by 568 nm line of Ar-Kr laser and by the 594 nm line of orange He-Ne laser. HiLyte Fluor™ 594 hydrazide is a carbonyl-reactive fluorescent labeling dye. It can be used for labeling glycoproteins such as HRP.
Spectra:	Maximum Ex/Em wavelength is 593/616 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-81271	HiLyte Fluor™ 594 acid
AS-81272-1 (1 mg)	HiLyte Fluor™ 594 acid, SE
AS-81272-5 (5 mg)	HiLyte Fluor™ 594 acid, SE

This product is for *in vitro* research use only