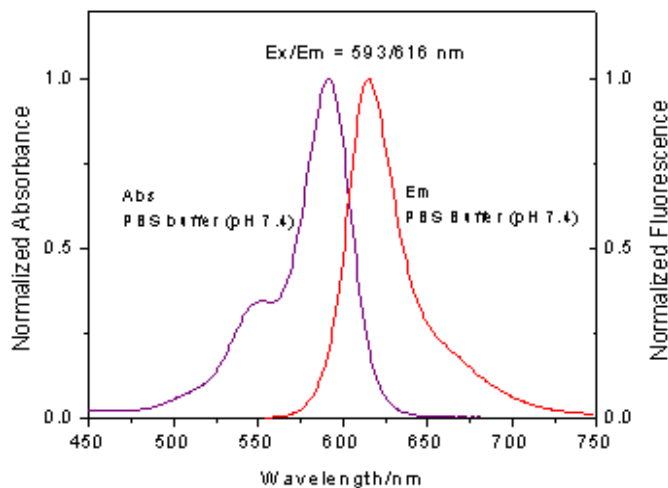




## Product Data Sheet

---

<b>Product Name:</b>	HiLyte Fluor™ 594 acid
<b>Size:</b>	10 mg
<b>Catalog Number:</b>	AS-81271
<b>Molecular Weight:</b>	751.87
<b>Solvent:</b>	DMF or DMSO
<b>Product Description:</b>	HiLyte Fluor™ 594, an alternative to Alexa Fluor™ 594 and DyLight Fluor™ 594, has spectral characteristics similar to those of Texas Red. The 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.
<b>Spectra:</b>	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-81272-1 (1 mg)	HiLyte Fluor™ 594 acid, SE
AS-81272-5 (5 mg)	HiLyte Fluor™ 594 acid, SE
AS-81274	HiLyte Fluor™ 594 hydrazide

This product is for *in vitro* research use only