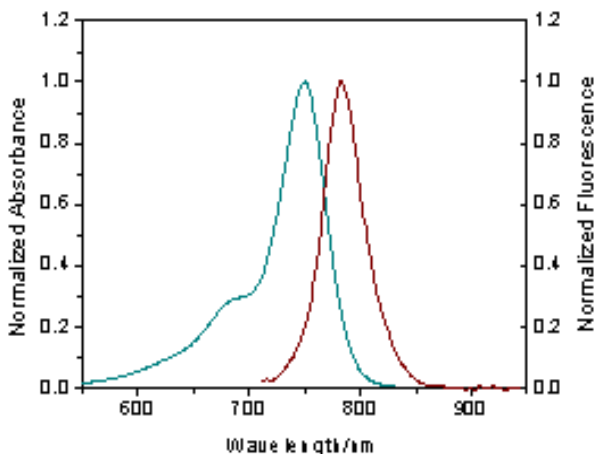




Product Data Sheet

Product Name:	HiLyte Fluor™ 750 C2 maleimide
Size:	1 mg
Catalog Number:	AS-81269
Molecular Weight:	1222.5
Solvent:	Water or DMF
Product Description:	Spectrally similar to Cy7 dye, HiLyte Fluor™ 750 C2 maleimide is the longest-wavelength thiol-reactive HiLyte Fluor™ dye currently available. Its fluorescence emission maximum at 778 nm is well separated from commonly used far-red fluorophores such as HiLyte Fluor™ 647, HiLyte Fluor™ 680 or allophycocyanin (APC), facilitating multicolor analysis. With a peak excitation at ~754 nm, conjugates of HiLyte Fluor™ 750 dyes are well excited by a xenon-arc lamp or dye-pumped lasers operating in the 720–750 nm range.
Spectra:	Maximum Ex/Em wavelength is 754/778 nm.



Stock Solution Preparation: Prepare 10 mg/ml dye stock solution in water or DMF. 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-81266	HiLyte Fluor™ 750 acid, SE
AS-81267	HiLyte Fluor™ 750 amine
AS-81268	HiLyte Fluor™ 750 hydrazide

This product is for *in vitro* research use only.