



## Product Data Sheet

---

**Product Name:** HiLyte Fluor™ 750 C2 maleimide

**Size:** 1 mg

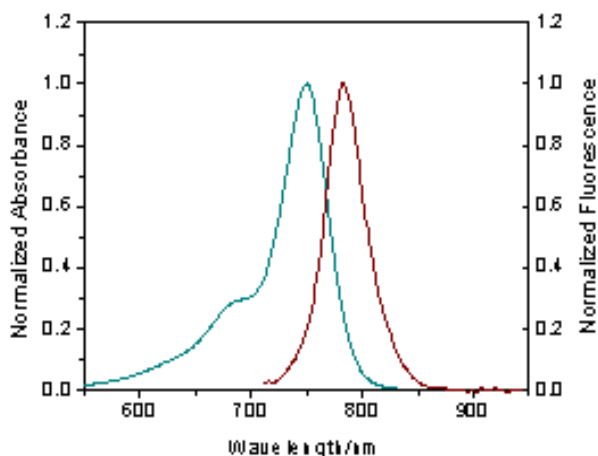
**Catalog Number:** 81269

**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
81266	HiLyte Fluor™ 750 acid, SE
81267	HiLyte Fluor™ 750 amine
81268	HiLyte Fluor™ 750 hydrazide

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