

## **Product Data Sheet**

Product Name: HiLyte Fluor™ 750 amine

Size: 1 mg

Catalog Number: OÈÜËB1267

Molecular Weight: 1071.38

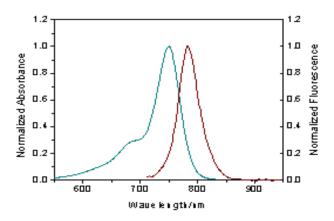
Solvent: Water

Product Description: HiLyte Fluor™ 750 amine is a carbonyl-reactive fluorescent labeling dye.

Spectrally similar to Cy7 dye, HiLyte Fluor<sup>TM</sup> 750 amine is the longest-wavelength carbonyl-reactive HiLyte Fluor<sup>TM</sup> dye currently available. Its fluorescence emission maximum at 778 nm is well separated from commonly used far-red fluorophores such as HiLyte Fluor<sup>TM</sup> 647, HiLyte Fluor<sup>TM</sup> 680 or Allophycocyanin (APC), facilitating multicolor analysis. With a peak excitation at ~754 nm, conjugates of HiLyte Fluor<sup>TM</sup> 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. 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
ŒÙËB1266	HiLyte Fluor™ 750 acid, SE
ŒÙËB126Ì	HiLyte Fluor™ 750 hydrazide
ŒÙËB1269	HiLyte Fluor™ 750 C2 maleimide

This product is for *in vitro* research use only