

Product Data Sheet

| Product Name: | HiLyte Fluor™ | 750 hydrazide |
|---------------|---------------|---------------|
|---------------|---------------|---------------|

Size: 1 mg

Catalog Number: ΆËB1268

Molecular Weight: 1043.32

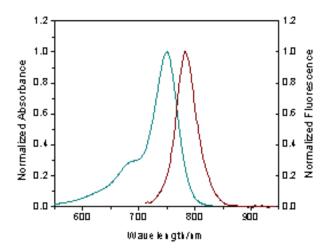
Solvent: Water or DMF

Product Description:

HiLyte Fluor[™] 750 hydrazide is a carbonyl-reactive fluorescent labeling dye. It can be used for labeling glycoproteins such as HRP. Spectrally similar to Cy7[™] dye, HiLyte Fluor[™] 750 hydrazide is the longest wavelength carbonyl-reactive HiLyte Fluor[™] dye currently available. Its fluorescence emission maximum at 782 nm is well separated from commonly used far-red fluorophores such as HiLyte Fluor[™] 647, HiLyte Fluor[™] 680 or Allophycocyanin (APC), facilitating multicolor analysis.

Spectra:

Maximum Ex/Em wavelength is 753/782 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 |
|-----------|--------------------------------|
| OEÙËB1266 | HiLyte Fluor™ 750 acid, SE |
| OEÙËB1267 | HiLyte Fluor™ 750 amine |
| OEÙËB1269 | HiLyte Fluor™ 750 C2 maleimide |

This product is for in vitro research use only