



## Product Data Sheet

---

**Product Name:** HiLyte™ Fluor 750 acid, SE

**Size:** 1 mg

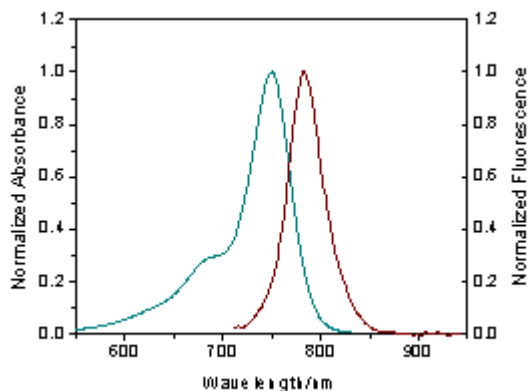
**Catalog Number:** AS-81266

**Molecular Weight:** 1328.75

**Solvent:** DMF or DMSO

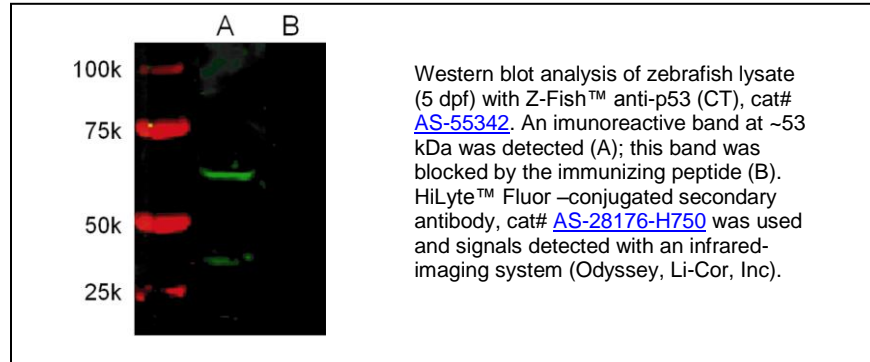
**Product Description:** Spectrally similar to Cy7™ dye, HiLyte™ Fluor 750 acid, SE dye is the longest-wavelength HiLyte™ Fluor dye currently available. Its fluorescence emission maximum at 778 nm is well separated from commonly used far-red fluorophores, including HiLyte™ Fluor 647, HiLyte™ Fluor 680 or allophycocyanin (APC), facilitating multicolor analysis. Extinction coefficient is 275,000 M<sup>-1</sup>cm<sup>-1</sup>, with a quantum yield of 0.12. With a peak excitation at ~753 nm, conjugates of HiLyte™ Fluor 750 dyes can be excited by a xenon-arc lamp or dye-pumped lasers operating in the 720–750 nm range. Its near infrared spectra render the HiLyte™ Fluor 750 dye an ideal candidate for *in vitro* and *in vivo* tissue imaging. HiLyte™ Fluor 750-protein conjugates are suitable for immunofluorescent staining, *in situ* hybridization, flow cytometry and other biological applications.

**Spectra:** Maximum Ex/Em wavelength is 754/778 nm.

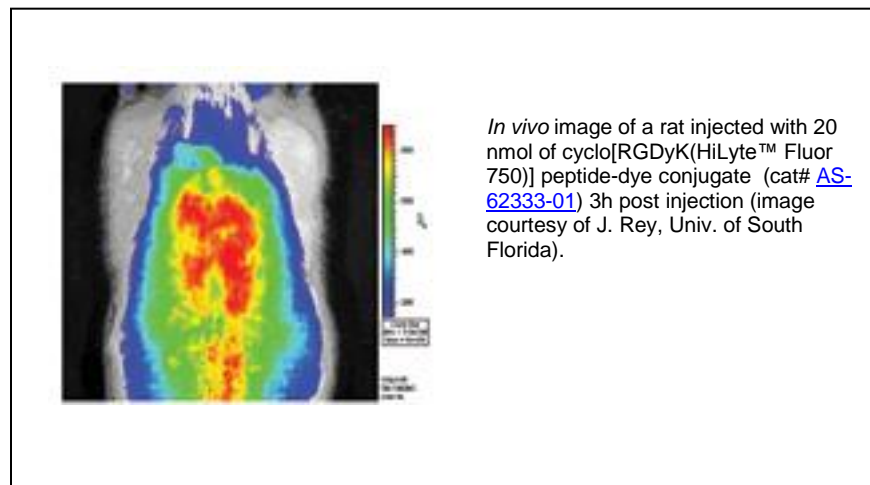


**Application:**

Western blot analysis with HiLyte™ Fluor 750 labeled secondary antibody



*In vivo* animal imaging with HiLyte™ Fluor 750 labeled RGD peptide



**Stock solution preparation:** Make 10 mg/ml of dye stock solution (7.5 mM) in high quality anhydrous DMSO. Completely dissolve all dye content by vortexing. Dye solution should be prepared fresh immediately before an experiment. Extensive storage of the dye solution 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-81267	<a href="#">HiLyte™ Fluor 750 amine</a>
AS-81268	<a href="#">HiLyte™ Fluor 750 hydrazide</a>
AS-81269	<a href="#">HiLyte™ Fluor 750 C2 maleimide</a>
AS-72043	<a href="#">AnaTag™ HiLyte™ Fluor 750 Protein Labeling Kit *Ultra Convenient</a>
AS-72044	<a href="#">AnaTag™ HiLyte™ Fluor 750 Microscale Protein Labeling Kit *Ultra Convenient</a>
	<a href="#">HiLyte™ Fluor 750 labeled Secondary Antibodies &amp; Streptavidin</a>

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