



Product Data Sheet

Product Name:	CyLyte Fluor 5, acid
Size:	5 mg, 25 mg, 100 mg
Catalog Number:	AS-89354-5, AS-89354-25, AS-89354-100
Molecular Weight:	656.81
CAS:	146368-11-8
Solvent:	DMF or water
Product Description:	CyLyte Fluor 5 dye has the same organic dye structure as that of Cy [®] 5. The CyLyte Fluor 5 dye is a far-red fluorescent label suitable for protein and nucleic acid conjugations used widely in imaging, immunocytochemistry, flow cytometry and FRET applications. It has significant advantages owing to its longer spectral wavelength over other fluorophores in minimizing the effect of potentially autofluorescent biological specimens and having pH insensitivity.
Spectral properties:	Abs/Em= 648/663 nm; $\epsilon = 250,000 \text{ M}^{-1}\text{cm}^{-1}$ in PBS buffer (pH7.4)
Stock Solution Preparation:	Prepare 10 mg/ml dye stock solution in DMF & 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 2 years if stored at -20°C, desiccated and protected from light.

This product is for *in vitro* research use only

Cy[®]5 is a registered trademark of GE Healthcare