



Product Data Sheet

Product Name:	CyLyte Fluor 3, acid
Size:	5 mg, 25 mg, 100 mg
Catalog Number:	AS-89353-5, AS-89353-25, AS-89353-100
Molecular Weight:	630.77
CAS:	146368-13-0
Solvent:	DMF or water
Product Description:	CyLyte Fluor 3 dye has the same organic dye structure as that of Cy [®] 3. The CyLyte Fluor 3 dye is a bright orange fluorescent dye suitable for protein and nucleic acid labeling used widely in imaging, immunocytochemistry, flow cytometry and FRET applications.
Spectral Properties:	Abs/Em= 550/564 nm; $\epsilon = 150,000 \text{ M}^{-1}\text{cm}^{-1}$ in PBS buffer (pH7.4)
Stock Solution Preparation:	Prepare 10 mg/ml dye stock solution in DMF or 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[®]3 is a registered trademark of GE Healthcare