



Product Data Sheet

| | |
|-----------------------------|---|
| Product Name: | CyLyte Fluor 5, NHS ester |
| Size: | 1 mg, 5 mg, 25 mg |
| Catalog Number: | AS-89357-1, AS-89357-5, AS-89357-25 |
| Molecular Weight: | 753.88 |
| CAS: | 146368-14-1 |
| Solvent: | DMF or DMSO |
| Product Description: | CyLyte Fluor 5 NHS ester dye has the same organic dye structure as that of Cy [®] 5 NHS ester dye. However, the counterion used in the final formulation of AnaSpec's CyLyte Fluor 5 NHS ester dye does vary compared to the Cy [®] 5 NHS ester dye. This CyLyte Fluor 5 NHS ester is a reactive dye used for the labeling of amino-groups in peptides, proteins, and oligonucleotides. This highly sensitive fluorophore has its emission maximum in the red region exhibiting maximum sensitivity thus minimizing background interference from biological specimen. |
| 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/DMSO. 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.