



Product Data Sheet

Product Name:	CyLyte Fluor 7, NHS ester
Size:	1 mg, 5 mg, 25 mg
Catalog Number:	AS-89358-1, AS-89358-5, AS-89358-25
Molecular Weight:	779.92
CAS:	477908-53-5
Solvent:	DMF or DMSO
Product Description:	CyLyte Fluor 7 NHS ester dye has the same organic dye structure as that of Cy [®] 7 NHS ester dye. However, the counterion used in the final formulation of AnaSpec's CyLyte Fluor 7 NHS ester dye does vary compared to the Cy [®] 7 NHS ester dye. The CyLyte Fluor 7 NHS ester is a near infrared amine reactive Cyanine 7 dye. This fluorophore can be used in conjugations for imaging applications in the near infrared region.
Spectral properties:	Abs/Em= 750/773 nm; $\epsilon = 270,000 \text{ M}^{-1}\text{cm}^{-1}$ in PBS buffer (pH7.4)
Stock Solution Preparation:	Prepare 10 mg/ml dye stock solution in DMF or 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[®]7 is a registered trademark of GE Healthcare.