

Product Data Sheet

Product Name: CyLyte Fluor 3, NHS ester

Size: 1 mg, 5 mg, 25 mg

Catalog Number: AS-89356-1, AS-89356-5, AS-89356-25

Molecular Weight: 727.85

CAS: 146368-16-3

Solvent: DMF or DMSO

Product Description: CyLyte Fluor 3 NHS ester dye has the same organic dye

structure as that of Cy®3 NHS ester dye. However, the counterion used in the final formulation of AnaSpec's CyLyte Fluor 3 NHS ester dye does vary compared to the Cy®3 NHS ester dye. The CyLyte Fluor 3 NHS ester is a reactive dye suitable for the labeling of amino-groups in biomolecules including soluble

proteins, peptides, and oligonucleotides/DNA.

Spectral Properties: Abs/Em= 550/564 nm; $\varepsilon = 150,000 \text{ M}^{-1}\text{cm}^{-1}$ in PBS buffer (pH7.4)

Stock Solution Preparation: Prepare 10 mg/ml dye stock solution in DMF or 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®3 is a registered trademark of GE Healthcare.