

Product Information Sheet

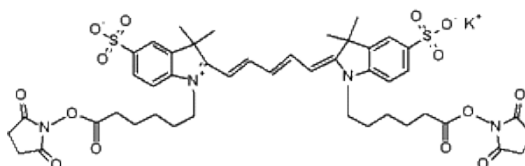
Ordering Information

Product Number:	282
Product Name:	Cyanine 5 bisuccinimidyl ester, potassium salt [same as GE Cy5 [®] bisNHS ester]
Unit Size:	1 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 975.13

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 649

Emission Wavelength: 665

Application Notes

This Cy5[®] dye is the same molecule to GE's Cy5[®] bisNHS ester. It readily reacts with amino groups. It has two NHS groups that might give brighter signal on some proteins than the Cy5 mono NHS ester (#151 and #280). Our Cy5[®] Fluors are thoroughly QC tested to ensure high levels of chromophore and reactive dye content. A variety of cyanine 5 (Cy5[®]) dyes has been used to label biological molecules for fluorescence imaging and other fluorescence-based biochemical analysis. They are widely used for labeling peptides, proteins and oligos etc. Cy5[®] dyes are one type of the most common red fluorophores. These versatile fluorophores can tolerate a pH range of 3-10 for use in a variety of applications at biologically relevant pHs. The dyes are also DMSO tolerant and photostable to enable transfer from storage to assay without loss of performance. The aqueous solubility eliminates the need for organic solvents in the assay buffers. Cy5[®] is the trademark of GE Healthcare.