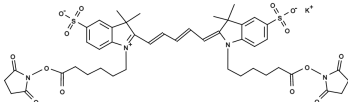


**Cyanine 5 bissuccinimidyl ester, potassium salt [same as GE Cy5® bisNHS ester]**Catalog number: 282  
Unit size: 1 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

**Chemical Properties**

Appearance	Dark blue solid
Molecular Weight	975.13
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	651 nm
Emission Wavelength	670 nm

**Applications**

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.