

## Cyanine 5 monosuccinimidyl ester, potassium salt [same as GE Cy5<sup>®</sup> NHS ester]

Catalog number: 280 Unit size: 1 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Blue solid
Molecular Weight	791.97
Soluble In	DMSO
Chemical Structure	$^{\circ}O_{3}S$ $\rightarrow$ $^{\circ}V$ $^{\circ$
Spectral Properties	
Excitation Wavelength	651 nm

## Applications

**Emission Wavelength** 

This Cy5<sup>®</sup> dye is the same molecule to GE's monoreactive Cy5<sup>®</sup> NHS ester. It readily reacts with amino groups. Our Cy5<sup>®</sup> Fluors are thoroughly QC tested to ensure high levels of chromophore and reactive dye content. Mono-reactive dyes are suitable for targeted, precise labeling of proteins and oligonucleotides and bis-reactive dyes are more suitable for general labeling. NHS ester dyes are recommended for labeling amine groups and maleimide dyes are recommended for labeling thiol groups. A variety of cyanine 5 (Cy5<sup>®</sup>) 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<sup>®</sup> 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<sup>®</sup> is the trademark of GE Healthcare.

670 nm