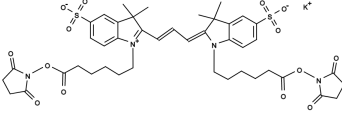


**Cyanine 3 bissuccinimidyl ester, potassium salt [same as GE Cy3® bisNHS ester]**Catalog number: 270  
Unit size: 1 mg**Product Details**

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

**Chemical Properties**

Appearance	Solid
Molecular Weight	949.10
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	555 nm
Emission Wavelength	569 nm

**Applications**

This Cy3® dye is the same molecule to GE's Cy3® bisNHS ester. It has two NHS groups that might give brighter signal on some proteins than the Cy3 mono NHS ester (#141 and #271). A variety of cyanine 3 (Cy3®) 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. Cy3® dyes have enhanced fluorescence upon binding to proteins. Cy3® NHS ester readily reacts with amino groups. Cy3® is the trademark of GE Healthcare.