

Product Information Sheet

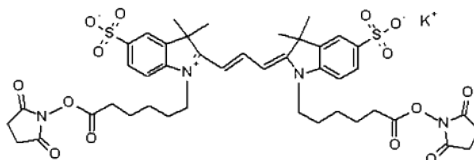
Ordering Information

| | |
|---------------------|---|
| Product Number: | 270 |
| Product Name: | Cyanine 3 bisuccinimidyl ester, potassium salt [same as GE Cy3 [®] 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: 949.10

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 555

Emission Wavelength: 565

Application Notes

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.