

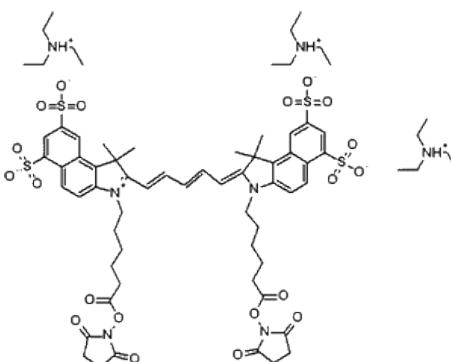
Product Information Sheet

Ordering Information

Product Number: 158
Product Name: Cyanine 5.5, bissuccinimidyl ester
Unit Size: 1 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 1500.87
Appearance: Blue solid
Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 678 nm
Emission Wavelength: 701 nm

Application Notes

A variety of cyanine 5.5 (Cy5.5®) 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. Cy5.5® dyes are one type of the most common red fluorophores. Cy5.5® NHS ester readily reacts with amino groups. Cy5.5® is the trademark of GE Healthcare.