

Product Information Sheet

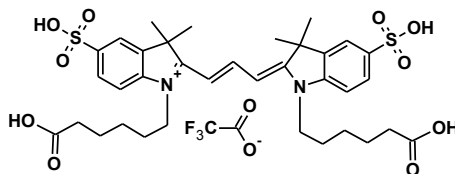
Ordering Information

Product Number:	137
Product Name:	Cyanine 3, bisacid
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight:	830.88
Appearance:	Purple solid

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	555 nm
Emission Wavelength:	565 nm

Application Notes

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.