

Product Information Sheet

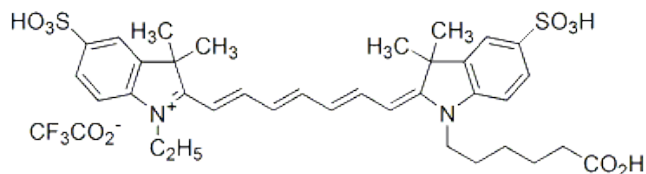
Ordering Information

Product Number:	160
Product Name:	Cyanine 7 monoacid [equivalent to Cy7 [®] acid]
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 796.87

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	749 nm
Emission Wavelength:	776 nm

Application Notes

A variety of cyanine 7 (Cy7[®]) 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. Cy7[®] dye conjugates are one type of the most common near infrared red fluorophores used in *in vivo* imaging applications. Cy7[®] is the trademark of GE Healthcare.