

Product Information Sheet

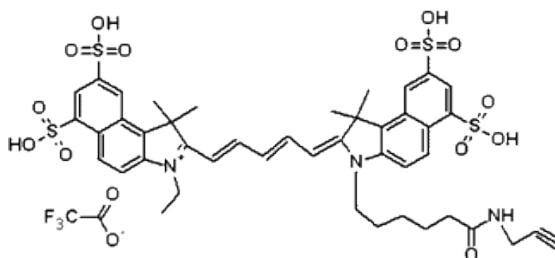
Ordering Information

Product Number:	179
Product Name:	Cyanine 5.5 alkyne [equivalent to Cy5.5 [®] alkyne]
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: 1068.14

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[®] alkyne selectively reacts with azide groups. Cy5.5[®] is the trademark of GE Healthcare.