

## Product Information Sheet

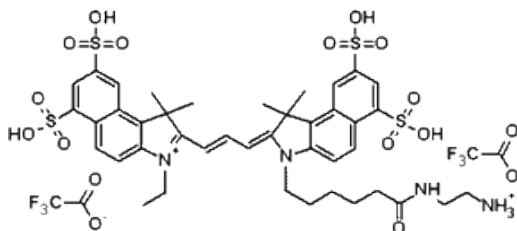
### Ordering Information

Product Number:	139
Product Name:	Cyanine 3.5 amine [equivalent to Cy3.5 <sup>®</sup> amine]
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: 1161.15

Chemical Structure:



Soluble in:	DMSO
Excitation Wavelength:	581 nm
Emission Wavelength:	596 nm

### Application Notes

A variety of cyanine 3.5 (Cy3.5<sup>®</sup>) 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. Cy3.5<sup>®</sup> dyes have enhanced fluorescence upon binding to proteins. Cy3.5<sup>®</sup> NHS ester readily reacts with carbonyl groups. Cy3.5<sup>®</sup> is the trademark of GE Healthcare.