

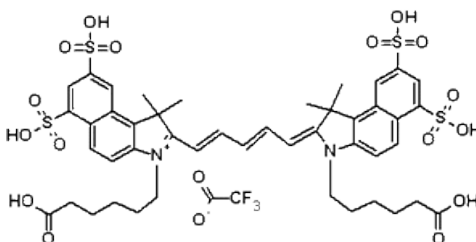
## Product Information Sheet

### Ordering Information

Product Number:	172
Product Name:	Cyanine 5.5 bis acid [equivalent to Cy5.5 <sup>®</sup> 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:	1117.17
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<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. Cy5.5<sup>®</sup> dyes are one type of the most common red fluorophores. Cy5.5<sup>®</sup> is the trademark of GE Healthcare.