

## Cyanine 5.5 bisacid [equivalent to Cy5.5® bisacid]

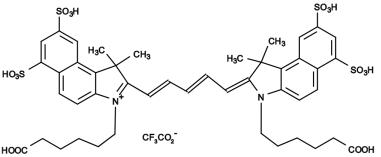
Catalog number: 172

Unit size: 5 mg

### Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

Appearance	Blue solid
Molecular Weight	1117.17
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

Excitation Wavelength	683 nm
Emission Wavelength	703 nm

### Applications

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® is the trademark of GE Healthcare.