

# Cyanine 7 monoacid [equivalent to Cy7® acid]

Catalog number: 160

Unit size: 5 mg

### **Product Details**

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

### **Chemical Properties**

Appearance Green solid

Molecular Weight 796.87

Soluble In DMSO

**Chemical Structure** 

### **Spectral Properties**

Excitation Wavelength 756 nm

Emission Wavelength 779 nm

## **Applications**

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.