

**Cy3B DBCO**Catalog number: 945  
Unit size: 1 mg**Product Details**

---

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

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	1154.39
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	560 nm
Emission Wavelength	571 nm

**Applications**

---

Cy3® dye is one of the most common cyanine dyes used for labeling proteins, nucleic acids and other biological molecules. A variety of Cy3® 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. However, Cy3 suffers low fluorescence intensity. It is the least fluorescent dyes among all the Cy dyes. Cy3B is an improved version of Cy3 dyes with significantly increased fluorescence quantum yield and photostability. Cy3B DBCO can be used for conjugations with azide-containing biological molecules under copper-free conditions. Cy3® is the trademark of GE Healthcare.