

## iFluor® IRD 800 CW SE

Catalog Number: 71610, 71611

Unit Size: 1 mg, 10 mg

### Product Details

---

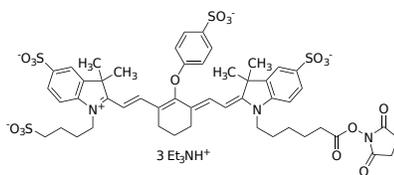
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid green
Molecular Weight	1403.83
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	774 nm
Emission Wavelength	792 nm

### Applications

---

iFluor® IRD 800 CW SE has the same chemical structure as the IRDye® 800 CW SE (IRDye® is the trademark of LI-COR). IRD 800CW SE is an amine reactive near-infrared (NIR) fluorescent dye with good water solubility. It can be used to label peptides, antibodies, and other amino-containing biomolecules such as amino-modified oligos. The IRD 800CW conjugates can be used as a probe for Western blot, flow cytometry, immunocytochemical assays, including In-Cell Western assays and On-Cell Western cell-based assays, protein arrays, microscopy, tissue section imaging, and *in vivo* imaging. IRD 800CW dye-conjugated probes are also used for fluorescence imaging-based cancer diagnosis.