

iFluor® IRD 800 CW Maleimide

Catalog Number: 71612 71613,

Unit Size: 1 mg 5 mg,

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	1239.33
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

iFluor® IRD 800 CW Maleimide has the same chemical structure as the IRDye® 800 CW Maleimide (IRDye® is the trademark of LI-COR). IRD 800CW Maleimide is a thiol-reactive near-infrared (NIR) fluorescent dye with good water solubility. It can be used to label peptides, antibodies, and other thiol-containing biomolecules such as thiol-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.