

## IVISense™ 680 NHS Fluorescent Dye (1mg)

Product Number: NEV11119

### DESCRIPTION

IVISense™ 680 NHS Fluorescent Dye is an amine reactive near-infrared (NIR) fluorochrome (NHS ester) for labeling biomolecules for in vivo imaging applications.

### MATERIAL

Available as a solid in 1 mg or 5 mg vials.

### STORAGE & HANDLING

- Upon receipt, IVISense™ 680 NHS Fluorescent Dye should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ 680 NHS Fluorescent Dye is stable for up to 12 months from the date of shipment.
- Allow to equilibrate to room temperature before opening.

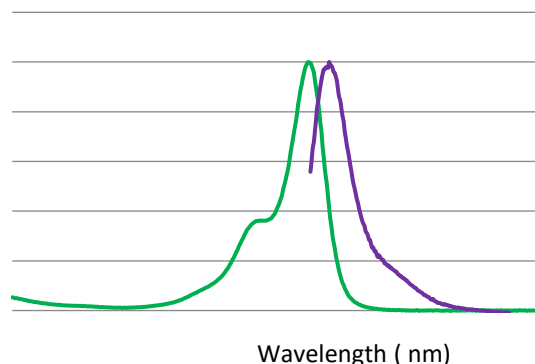
### IN VIVO IMAGING

- Optimized in vivo NIR label for labeling target molecules including peptides, small molecules, nanoparticles, and macromolecules.
- Highly hydrophilic and suitable for single or multiple dye loading.

Property	Specification
MW	1856 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	665 ± 5 nm
• Emission	688 ± 5 nm
Absorbance <sup>1</sup>	668 ± 3 nm
Extinction	~210,000 M <sup>-1</sup> cm <sup>-1</sup>
Purity <sup>2</sup>	>95%
Appearance	Blue Solid

1. Absorbance and fluorescence maxima in 1xPBS.
2. Reactive ester as determined by RP-HPLC, measuring absorbance at 670 nm

\*MW as Penta TEA Salt



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS, OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

NEV11119-R Rev01