

IVISense™ 800 NHS Fluorescent Dye (1mg)

Product Number: NEV11107

DESCRIPTION

IVISense™ 800 NHS Fluorescent Dye is an amine reactive (NHS Ester) near-infrared (NIR) fluorochrome 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™ 800 NHS Fluorescent Dye should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ 800 NHS Fluorescent Dye is stable for up to 12 months from the date of shipment.
- Allow to equilibrate to room temperature before opening.
- Recommended solvents for reconstitution:
Aqueous buffers pH 6.0-7.0
Organic solvents: DMSO, DMF
- Recommended solvents for performing conjugation reactions:
Aqueous buffers pH 7.5-8.5
Organic solvents: DMSO, DMF
Carbonate buffers are not recommended.

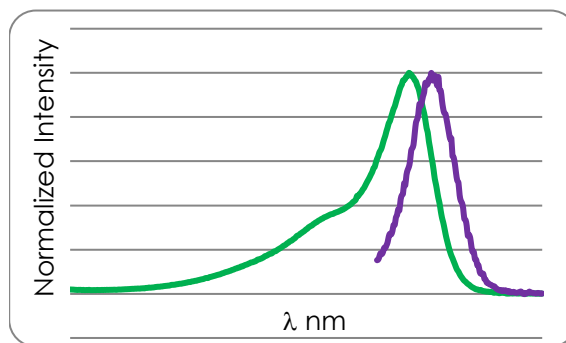
IN VIVO IMAGING

- Optimized in vivo NIR label for labeling target molecules including peptides, small molecules, macromolecules, and nanoparticles.

Property	Specification
MW	1464 g mol ⁻¹
Fluorescence ¹	
• Excitation	785 nm
• Emission	810 nm
Absorbance ¹	787 ± 5 nm
Extinction	~200,000 M ⁻¹ cm ⁻¹
Purity ²	>90%
Appearance	Green Solid

1. Absorbance and fluorescence maxima in 1xPBS.
2. Reactive ester as determined by RP-HPLC, measuring absorbance at 790 nm

*MW as Tris 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
NEV11107-R Rev01