

IVISense™ 680 NHS Fluorescent Self-Quenching Dye (1mg)

Product Number: NEV10121

DESCRIPTION

IVISense™ 680 NHS Fluorescent Self-Quenching 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 Self-Quenching Dye should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ 680 NHS Fluorescent Self-Quenching Dye is stable for up to 24 months from the date of shipment.
- Allow to equilibrate to room temperature before opening.

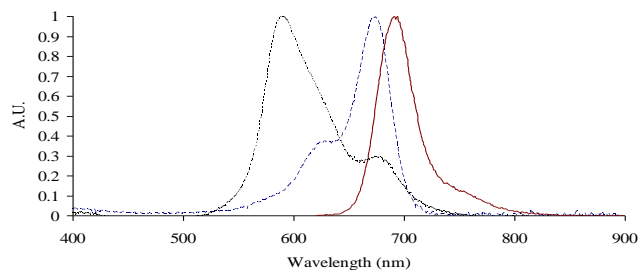
IN VIVO IMAGING

- Optimized in vivo NIR label for labeling target molecules including peptides, small molecules, macromolecules, and nanoparticles.
- Labeling of target molecules that have multiple amine reactive sites with > 1 fluorochrome results in a blue shift in the absorbance spectra.

Property	Specification
MW	1439 g mol ⁻¹
Fluorescence ¹	
• Excitation	673 ± 5 nm
• Emission	691 ± 5 nm
Absorbance ¹	673 ± 5 nm
Extinction	~230,000 M ⁻¹ cm ⁻¹
Purity ²	>90%
Appearance	Blue Solid

1. Absorbance and fluorescence maxima in 1xPBS.
2. As determined by RP-HPLC, measuring absorbance at 675nm

*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
NEV10121-R Rev01