

IVISense™ Acute Vascular 680 Fluorescent Probe (5mg)

Product Number: NEV10130

DESCRIPTION

IVISense™ Acute Vascular 680 Fluorescent Probe is the acid form of IVISense™ 680 Fluorescent Dye that can be used as a control or a fluorescent in vivo blood pool imaging agent. Upon dilution, a 5mg package of IVISense™ Acute Vascular 680 Fluorescent Probe provides sufficient reagents for imaging approximately 20 mice (weighing ~25 grams each) when using the recommended dose of 0.05mg/100 μ L 1xPBS.

MATERIAL

Available as a solid in 5 mg vials.

STORAGE & HANDLING

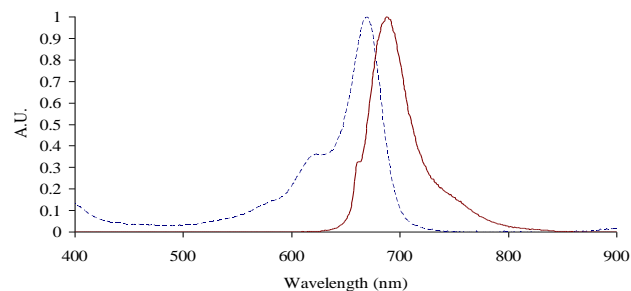
- Upon receipt, IVISense™ Acute Vascular 680 Fluorescent Probe should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Acute Vascular 680 Fluorescent Probe is stable for up to 24 months from the date of shipment.
- Allow to equilibrate to room temperature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

IN VIVO IMAGING

- The recommended procedure for in vivo imaging with IVISense™ Acute Vascular 680 Fluorescent Probe is administration via tail vein injection and imaging 30 minutes post injection.
- Imaging in materigel plugs and tumor models: IVISense™ Tomato Lectin 680 Fluorescent Probe can be used as a control during Imaging studies as a blood pool imaging agent.

Property	Specification
MW	~1646 g mol ⁻¹
Fluorescence ¹	
• Excitation	670 ± 5 nm
• Emission	688 ± 5 nm
Absorbance ¹	670 ± 5 nm
Extinction	~210,000 M ⁻¹ cm ⁻¹
Purity ²	>95%
Appearance	Blue Solid

1. Absorbance and fluorescence maxima in 1xPBS.
 2. As determined by RP-HPLC, measuring absorbance at 670nm
- *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
NEV10130-R Rev01