

IVISense™ Acute Vascular 750 Fluorescent Probe (5mg)

Product Number: NEV10177

DESCRIPTION

IVISense™ Acute Vascular 750 Fluorescent Probe is the acid form of IVISense™ 750 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 750 Fluorescent Probe provides sufficient reagents for imaging approximately 20 mice (weighing ~25 grams each) when using the recommended dose of 0.05mg/100 µL 1xPBS.

MATERIAL

Available as a solid in 5 mg vials.

STORAGE & HANDLING

- Upon receipt, IVISense™ Acute Vascular 750 Fluorescent Probe should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Acute Vascular 750 Fluorescent Probe is stable for up to 24 months from the date of shipment.
- Allow to equilibrate to room temperature before opening.

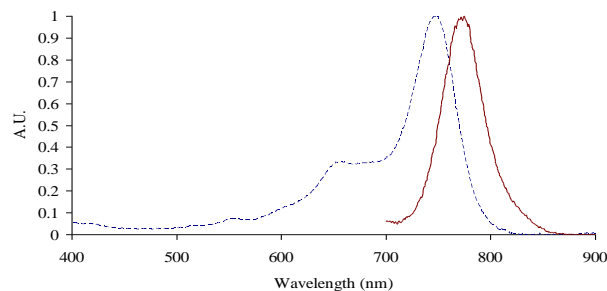
IN VIVO IMAGING

- Dilute IVISense™ Acute Vascular 750 to a concentration of 0.5mg/mL with sterile 1xPBS.
- The recommended procedure for in vivo imaging with IVISense™ Acute Vascular 750 Fluorescent Probe is administration via tail vein injection and imaging 30 minutes post injection.
- IVISense™ Acute Vascular 750 can be used as a control during imaging studies and as a blood pool imaging agent.

Property	Specification
MW (acid form)	1086 g mol ⁻¹
Fluorescence ¹	
• Excitation	750 ± 5 nm
• Emission	775 ± 5 nm
Absorbance ¹	750 ± 5 nm
Extinction	~240,000 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 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
NEV10177-R Rev01