

IVISense™ Edema 680 Fluorescent Probe

Product Number: NEV10116

DESCRIPTION

IVISense™ Edema 680 Fluorescent Probe is a small molecule fluorescent in vivo blood pool imaging agent. IVISense™ Edema 680 Fluorescent Probe enables imaging of circulation, blood vessels, vasculature, vascular leak, including that associated with early oncologic and ophthalmologic lesions.

MATERIAL

Each vial contains 40 nmoles of IVISense™ Edema 680 Fluorescent Probe in 1500 ul at a concentration of 4 nmol/150 μL, in 1xPBS. The IVISense™ Edema 680 Fluorescent Probe solution has been filtered through a 0.2 μm filter. This material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 4 nmoles of IVISense™ Edema 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

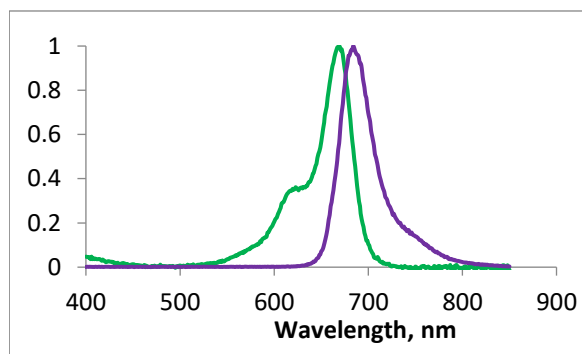
- Upon receipt, IVISense™ Edema 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Edema 680 Fluorescent Probe is stable for up to 24 months from the date of shipment.
- Allow IVISense™ Edema 680 Fluorescent Probe imaging agent to equilibrate to room temperature before injecting into animals.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Edema 680 Fluorescent Probe is administration via tail vein injection and imaging 0.5-24 hours post injection.
- IVISense™ Edema 680 Fluorescent Probe enables imaging of circulation, blood vessels, vasculature, vascular leak, including that associated with early oncologic and ophthalmologic lesions.

Property	Specification
MW	~1540 g mol ⁻¹
Fluorescence ¹	
• Excitation	675 ± 5 nm
• Emission	692 ± 5 nm
Absorbance ¹	675 ± 5 nm
Purity ²	>95%
Appearance	Clear blue solution

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



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
NEV10116-R Rev01