

IVISense™ Vascular 680 Fluorescent Probe

Product Number: NEV10054EX

DESCRIPTION

IVISense™ Vascular 680 Fluorescent Probe is a fluorescent in vivo blood pool imaging agent. IVISense™ Vascular 680 Fluorescent Probe is a near-infrared labeled fluorescent macromolecule that remains localized in the vasculature for extended periods of time and enables imaging of blood vessels and angiogenesis..

MATERIAL

Each vial contains 24 nmol of IVISense™ Vascular 680 Fluorescent Probe as a lyophilized solid. The solution of IVISense™ Vascular 680 Fluorescent Probe has been filtered through a 0.2 µm filter prior to lyophilization. Upon reconstitution with 1.2 mL of 1 x PBS, this material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/100 µL 1xPBS of IVISense™ Vascular 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

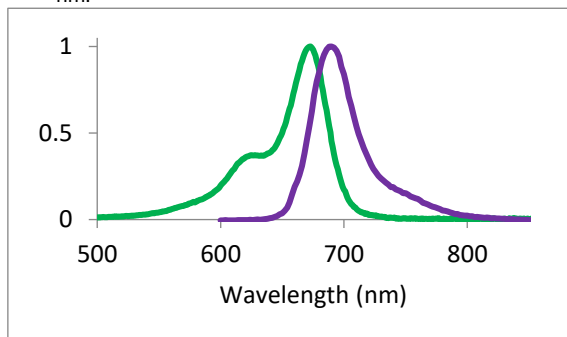
- Upon receipt, IVISense™ Vascular 680 Fluorescent Probe should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Vascular 680 Fluorescent Probe is stable for up to six months from the date of shipment.
- Once reconstituted, the PBS solution is stable up to 14 days when stored at 2-8 °C.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Vascular 680 Fluorescent Probe is administration via tail vein injection and imaging 24 hours post injection.
- Imaging in Arthritis: IVISense™ Vascular 680 Fluorescent Probe can be used to characterize vascular changes and therapeutic responses associated with animal models of arthritis.
- Imaging in Oncology: IVISense™ Vascular 680 Fluorescent Probe can be used to study angiogenesis, as a marker for blood vessel density, in animal tumor models.

| Property | Specification |
|---------------------------|------------------------------|
| MW | ~ 70,000 g mol ⁻¹ |
| Fluorescence ¹ | |
| • Excitation | 670 nm |
| • Emission | 690 nm |
| Absorbance ¹ | 673 ±10nm |
| Purity ² | >90% |
| Appearance | Blue solid |

1. Absorbance, excitation, and fluorescence maxima of IVISense™ Vascular 680 Fluorescent Probe in 1xPBS.
2. As determined by SE-HPLC and measuring absorbance at 680 nm.



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
 NEV10054EX-R Rev01