

IVISense™ Vascular NP 750 Fluorescent Nanoparticles

Product Number: NEV10150

DESCRIPTION

IVISense™ Vascular NP is a highly fluorescent near infrared nanoparticle specifically designed for in vivo imaging. IVISense Vascular NP 750 contains an iron oxide core that is coated to specifically produce a functionalized biocompatible probe comprised of a pegylated fluorescent nanoparticle that remains localized in the vasculature for extended periods of time and enables imaging of blood vessels, angiogenesis, blood-brain-barrier compromise, and vascular cell fluid-phase pinocytotic function.

MATERIAL

Each vial contains 500 µL of IVISense™ Vascular NP 750 Fluorescent Nanoparticles in 1xPBS. The IVISense™ Vascular NP 680 Fluorescent Nanoparticles solution has been filtered through a 0.2 µm filter. This material provides sufficient reagent for imaging approximately 5 mice (weighing ~25 grams each) when using the recommended injection volume of 100 µL of IVISense™ Vascular NP 750 Fluorescent Nanoparticles per mouse.

STORAGE & HANDLING

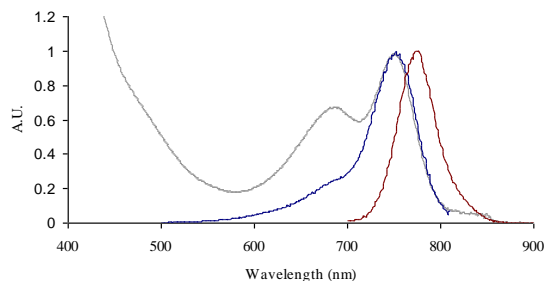
- Upon receipt, IVISense™ Vascular NP 750 Fluorescent Nanoparticles should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT. DO NOT FREEZE.
- When stored and handled properly, IVISense™ Vascular NP 750 Fluorescent Nanoparticles is stable for up to 12 months.
- Allow to equilibrate to room temperature before use.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for analysis by intravital microscopy imaging using IVISense™ Vascular NP 750 Fluorescent Nanoparticles is administration via tail vein injection and imaging 0 - 4 hour post injection
- IVISense™ Vascular NP 750 Fluorescent Nanoparticles can also be used at later time points (>12 hours) to image vascular leak in tissue site of inflammation and cancer. IVISense™ Vascular NP 680 Fluorescent Nanoparticles can be imaged within the interstitium for up to 24 hours post tail vein injection.

Property	Specification
Particle Size	20-50nm
Fluorescence ¹	
• Excitation	750 ± 5 nm
• Emission	775 ± 5 nm
Extinction ¹	4 X 10 ⁶ ± 2 X 10 ⁶ M ⁻¹ cm ⁻¹
Appearance	Greenish-brown solution

1. Absorbance, excitation, and fluorescence maxima of in 1xPBS.



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
NEV10150-R Rev01