

IVISense™ Integrin Receptor 750 Fluorescent Probe

Product Number: NEV10873

DESCRIPTION

IVISense™ Integrin Receptor 750 Fluorescent Probe is a targeted fluorescence imaging agent comprising a potent, selective non-peptide small molecule integrin $\alpha_v\beta_3$ antagonist and an NIR fluorochrome. This agent has been developed to enable in vivo visualization and quantification of integrin $\alpha_v\beta_3$ expression in neovasculature as well as in tumor cells, to monitor tumor angiogenesis, growth, and treatment efficacy. Half-life in tissue of IVISense™ Integrin Receptor 750 Fluorescent Probe signal is approximately 4 days.

MATERIAL

Each vial contains 24 nmol of IVISense™ Integrin Receptor 750 Fluorescent Probe lyophilized solid. The IVISense™ Integrin Receptor 750 Fluorescent Probe solution has been filtered through a 0.2 μm filter prior to lyophilization. Reconstitute IVISense™ Integrin Receptor 750 Fluorescent Probe with 1.2 mL of 1 x PBS before injecting into animals. The packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol (100 μL) of IVISense™ Integrin Receptor 750 Fluorescent Probe per mouse.

STORAGE & HANDLING

- Upon receipt, IVISense™ Integrin Receptor 750 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Integrin Receptor 750 Fluorescent Probe is stable for up to 12 months in the lyophilized form.
- Before opening the vial, ensure that the lyophilized powder is present at the bottom of the vial.
- Once reconstituted, the solution is stable up to 14 days when stored at 2-8°C and protected from light.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Integrin Receptor 750 Fluorescent Probe is administration via intravenous injection and imaging 24 hours post injection.
- Imaging maybe performed as early as 6 hours with some reduction in target signal/noise. It will clear from tissues after approximately 5. Repeat injection and imaging may be performed every 5 days for longitudinal studies.
- IVISense™ Integrin Receptor 750 Fluorescent Probe enables imaging of neovasculature and tumors in a range of oncology applications.

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
NEV10873-R Rev01

Property	Specification
MW	1278 g mol ⁻¹
Fluorescence ¹	
• Excitation	755 nm
• Emission	775 nm
Absorbance ¹	748 nm
Purity ²	>95%
Appearance	Dark blue-green solid

1. Absorbance, excitation, and fluorescence maxima of in 1xPBS.
2. As determined by RP-HPLC, measuring absorbance at 750nm

