

IVISense™ Integrin Receptor 680 Fluorescent Probe

Product Number: NEV10645

DESCRIPTION

IVISense™ Integrin Receptor 680 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 680 Fluorescent Probe signal is 24 hours.

MATERIAL

Each vial contains 24 nmol of IVISense™ Integrin Receptor 680 Fluorescent Probe lyophilized solid. The IVISense™ Integrin Receptor 680 Fluorescent Probe solution has been filtered through a 0.2 μm filter prior to lyophilization. Reconstitute IVISense™ Integrin Receptor 680 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 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

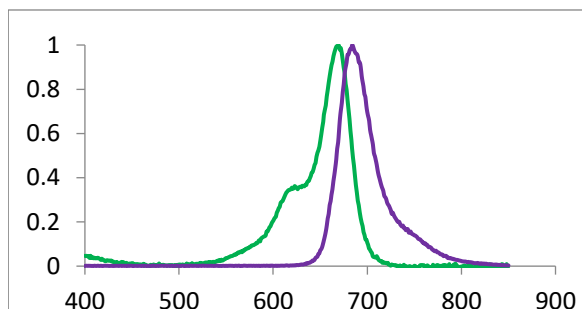
- Upon receipt, IVISense™ Integrin Receptor 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Integrin Receptor 680 Fluorescent Probe is stable for up to 12 months in the lyophilized form.
- Once reconstituted, the solution 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 680 Fluorescent Probe is administration via intravenous injection and imaging 3-48 hours post injection.
- Imaging at earlier time points (~3 hours) is recommended when imaging the vasculature.
- Imaging at later time points (24-48 hours) is recommended when imaging tumors to reduce background.
- IVISense™ Integrin Receptor 680 Fluorescent Probe enables imaging of neovasculature and tumors in a range of oncology applications.

Property	Specification
MW	1432 g mol ⁻¹
Fluorescence ¹	
• Excitation	675 nm
• Emission	693 nm
Absorbance ¹	675 ± 5 nm
Purity ²	>95%
Appearance	Dark blue solid

1. Absorbance, excitation, and fluorescence maxima of 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
NEV10645-R Rev01