

# IVISense™ Tomato Lectin 680 Fluorescent Probe

Product Number:      NEV10060

## DESCRIPTION

IVISense™ Tomato Lectin 680 Fluorescent Probe is a fluorescent in vivo endothelial cell imaging agent. IVISense™ Tomato Lectin 680 Fluorescent Probe is a near-infrared labeled fluorescent macromolecule that targets the vasculature and enables imaging of blood vessels and angiogenesis.

## MATERIAL

Each vial contains 24 nmol of IVISense™ Tomato Lectin 680 Fluorescent Probe as a lyophilized solid. The 1XPBS solution of IVISense™ Tomato Lectin 680 Fluorescent Probe has been filtered through a 0.2 µm filter prior to lyophilization. Upon reconstitution with 1.2 mL of DISTILLED WATER, this 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™ Tomato Lectin 680 Fluorescent Probe per mouse. The lyophilized formulation contains salt. When reconstituted with the recommended amount of water the salt concentration will be equivalent to 1XPBS.

## STORAGE & HANDLING

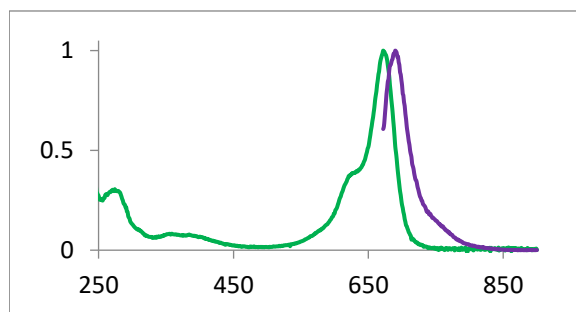
- Upon receipt, IVISense™ Tomato Lectin 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Tomato Lectin 680 Fluorescent Probe is stable for 3 months from the date of shipment.
- Once reconstituted, the solution is stable for up to 14 days when store at 2-8 °C and protected from light.

## IN VIVO IMAGING

- The recommended procedure for in vivo imaging with IVISense™ Tomato Lectin 680 Fluorescent Probe is administration via intravenous injection and imaging 6 hours post injection.
- Imaging in matrigel plugs and tumor models: IVISense™ Tomato Lectin 680 Fluorescent Probe can be used to study angiogenesis and blood vessel density, in matrigel plugs and animal tumor models.

Property	Specification
MW	~72,000 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	670 nm
• Emission	690 nm
Absorbance <sup>1</sup>	675 ± 10 nm
Purity <sup>2</sup>	>95%
Appearance	Blue Solid

1. Absorbance, excitation, and fluorescence maxima 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  
 NEV10060-R Rev01