

# IVISense™ Pan Cathepsin 750 Fluorescent Probe

Product Number:      NEV10001EX

## DESCRIPTION

IVISense™ Pan Cathepsin 750 Fluorescent Probe is a protease activatable fluorescent in vivo imaging agent that is activated by key disease associated proteases such as Cathepsin B, L, S and Plasmin. IVISense™ Pan Cathepsin 750 Fluorescent Probe is optically silent in its unactivated state and becomes highly fluorescent following protease-mediated activation.

## MATERIAL

Each vial contains 24 nmol (fluorophore) of IVISense™ Pan Cathepsin 750 Fluorescent Probe as a lyophilized solid. The solution of IVISense™ Pan Cathepsin 750 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™ Pan Cathepsin 750 Fluorescent Probe per mouse.

## STORAGE & HANDLING

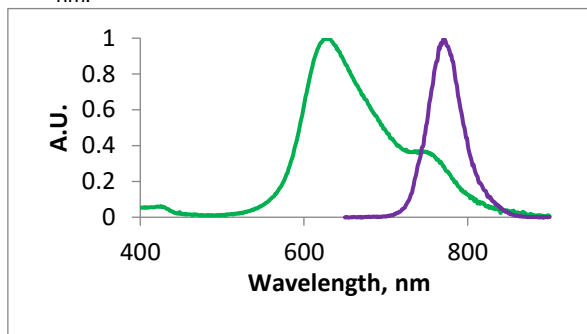
- Upon receipt, IVISense™ Pan Cathepsin 750 Fluorescent Probe should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Pan Cathepsin 750 Fluorescent Probe is stable for up to six months from the date of shipment.
- Once reconstituted, the PBS solution 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™ Pan Cathepsin 750 Fluorescent Probe is administration via tail vein injection and imaging 24 hours post injection.
- Imaging in Arthritis: IVISense™ Pan Cathepsin 750 Fluorescent Probe can be used as a marker for disease progression and therapeutic response in animal models of arthritis.
- Imaging in Oncology: IVISense™ Pan Cathepsin 750 Fluorescent Probe can be used as a marker for disease progression in animal tumor models.

Property	Specification
MW	~ 450,000 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	750 nm
• Emission	770 nm
Absorbance <sup>1</sup>	630 ±10nm
Purity <sup>2</sup>	>95%
Appearance	Blue solid

1. Absorbance, excitation, and fluorescence maxima of IVISense™ Pan Cathepsin 750 Fluorescent Probe in 1xPBS.
2. As determined by SE-HPLC and measuring absorbance at 750 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  
NEV10001EX-R Rev01