

IVISense™ Pan Cathepsin 680 Fluorescent Probe

Product Number: NEV10003

DESCRIPTION

IVISense™ Pan Cathepsin 680 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 680 Fluorescent Probe is optically silent in its unactivated state and becomes highly fluorescent following protease-mediated activation.

MATERIAL

Each vial contains 20 nmol of IVISense™ Pan Cathepsin 680 Fluorescent Probe at a concentration of 20 nmol/150 µL, in 1xPBS. The IVISense™ Pan Cathepsin 680 Fluorescent Probe solution has been filtered through a 0.2 µm filter. Upon dilution with 1350 µl of 1 x PBS, this material provides sufficient reagents for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/150 µL 1xPBS of IVISense™ Pan Cathepsin 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

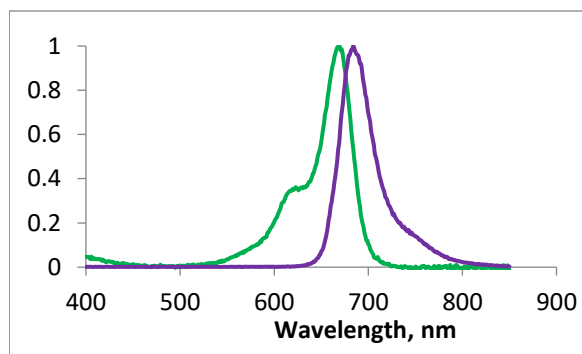
- Upon receipt, IVISense™ Pan Cathepsin 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Pan Cathepsin 680 Fluorescent Probe is stable for up to twelve months from the date of shipment.
- Allow IVISense™ Pan Cathepsin 680 Fluorescent Probe imaging agent to equilibrate to room temperature before injecting into animals.

IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Pan Cathepsin 680 Fluorescent Probe is administration via tail vein injection and imaging 24 hours post injection.
- Imaging in Arthritis: IVISense™ Pan Cathepsin 680 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 680 Fluorescent Probe can be used as a marker for disease progression in animal tumor model

Property	Specification
MW	~400,000 g mol ⁻¹
Fluorescence ¹	
• Excitation	680 ± 10 nm
• Emission	700 ± 10 nm
Absorbance ¹	680 ± 10 nm
Purity ²	>95%
Appearance	Clear blue solution

1. Absorbance, excitation, and fluorescence maxima of IVISense™ Pan Cathepsin 680 Fluorescent Probe 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
NEV10003-R Rev01