

IVISense™ MMP 680 Fluorescent Probe

Product Number: NEV10126

DESCRIPTION

IVISense™ MMP 680 Fluorescent Probe is a protease activatable fluorescent in vivo imaging agent that is activated by key matrix metalloproteinases (MMPs) including MMP-2, -3, -9 and -13. IVISense™ MMP 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™ MMP 680 Fluorescent Probe at a concentration of 20 nmol/1.5 mL in 1xPBS. The IVISense™ MMP 680 Fluorescent Probe solution has been filtered through a 0.2 µm filter. This package 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™ MMP 680 Fluorescent Probe per mouse.

STORAGE & HANDLING

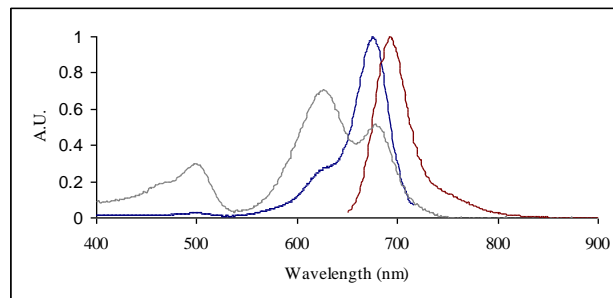
- Upon receipt, IVISense™ MMP 680 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ MMP 680 Fluorescent Probe is stable for up to six months from the date of shipment.
- Allow IVISense™ Pan Cathepsin 680 Fluorescent Probe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

IN VIVO IMAGING & APPLICATIONS

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

Property	Specification
MW	~450,000 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™ MMP 680 Fluorescent Probe in 1xPBS.
2. As determined by SE-HPLC, measuring absorbance at 680nm



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
NEV10126-R Rev01