

IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe

Product Number: NEV11169

DESCRIPTION

IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe is a member of a family of activatable fluorescent imaging agents comprising a novel architecture, termed F.A.S.T. (Fluorescent Activatable Sensor Technology), that confers an improved pharmacokinetic profile with earlier imaging time points. This architecture offers higher target specific signal with reduced background while also reducing the optimal imaging time after injection.

IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe is a neutrophil elastase activatable agent that is optically silent upon injection and produces fluorescent signal after cleavage by elastase produced by neutrophil cells.

Neutrophil elastase is a key protease involved in acute lung injury, acute respiratory distress syndrome, as well as many other inflammatory processes such as emphysema, cystic fibrosis, COPD, wound healing, rheumatoid arthritis, ischemia-reperfusion, and many others. IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe is a selective neutrophil elastase-activatable agent designed for imaging of this serine protease both in vitro and in vivo. The half-life in plasma of IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe is 4 hours.

MATERIAL

IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe: Each vial contains 48 nmol of IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe in dry solid form.

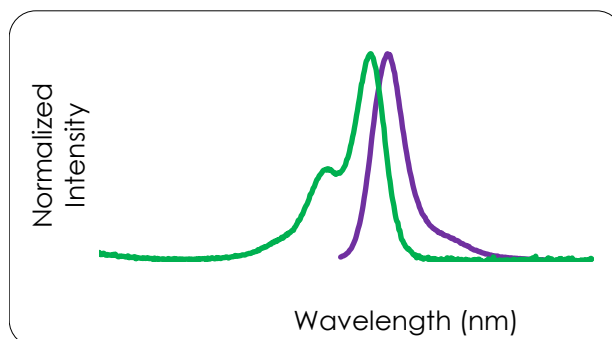
IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe has been filtered through a 0.2 µm filter prior to drying. Reconstitute IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe with 1.2 mL of 1 x PBS before injecting into animals. The packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 4 nmols (100 µL) of IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe per mouse.

STORAGE & HANDLING

- Upon receipt, IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense Neutrophil Elastase 680 FAST Fluorescent Probe is stable for up to 12 months in dry solid form.
- Before opening the vial, check to ensure that all of the solid material is at the bottom of the vial.

Property	Specification
MW	~43,000 g mol ⁻¹
Fluorescence ¹	
• Excitation	675 nm
• Emission	693 nm
Purity ²	>95%
Appearance	Blue Solid

1. Absorbance and fluorescence in 1xPBS.
2. As determined by RP-HPLC, measuring absorbance at 675nm



- 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 generally recommended procedure for in vivo imaging with IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe is administration via intravenous injection and imaging 4-8 hours post injection.
- IVISense™ Neutrophil Elastase 680 FAST Fluorescent Probe enables imaging of neutrophil elastase activity in applications including acute lung injury models, acute respiratory distress syndrome, emphysema, cystic fibrosis, COPD, wound healing, rheumatoid arthritis, and ischemia-reperfusion.

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
NEV11169-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.