

# IVISense™ Annexin-V 750 Fluorescent Probe

Product Number: NEV11053

## DESCRIPTION

IVISense™ Annexin-V 750 Fluorescent Probe is a targeted fluorescence imaging agent comprising a selective protein, Annexin A5 and a near-infrared (NIR) fluorochrome. This agent has been developed to enable in vivo visualization and quantification of the membrane-bound phospholipid, phosphatidylserine, exposed in the outer leaflet of the cell membrane lipid bilayer during the early stages of apoptosis.

## MATERIAL

Each vial contains 10 mouse doses of IVISense™ Annexin-V 750 Fluorescent Probe in solution. IVISense™ Annexin-V 750 Fluorescent Probe is supplied in 1000 µL of 25 mM HEPES/NaOH, 140 mM NaCl, pH 7.4. The packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 100 µL of IVISense™ Annexin-V 750 Fluorescent Probe per mouse.

## STORAGE & HANDLING

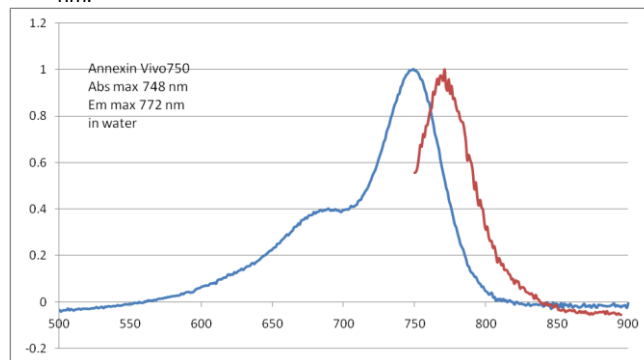
- Upon receipt, IVISense™ Annexin-V 750 Fluorescent Probe should be IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, IVISense™ Annexin-V 750 Fluorescent Probe is stable for six months from the date of shipment.

## IN VIVO IMAGING & APPLICATIONS

- The recommended procedure for in vivo imaging with IVISense™ Annexin-V 750 Fluorescent Probe is administration via intravenous injection and imaging 2 hours post injection.
- IVISense™ Annexin-V 750 Fluorescent Probe will clear from tissues after ~3 days after which repeat injection and imaging may be performed for longitudinal studies. However, it is recommended that a pre-injection baseline image be taken prior to reinjection and imaging.
- IVISense™ Annexin-V 750 Fluorescent Probe enables imaging of therapeutic response in oncology models, and has applications in cardiovascular disease, transplant rejection, and pancreatitis models.

Property	Specification
MW	35,000 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	755 nm
• Emission	772 nm
Absorbance <sup>1</sup>	748 nm
Purity <sup>2</sup>	>95%
Appearance	Blue solution

1. Absorbance and fluorescence maxima in water
2. As determined by RP-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  
NEV11053-R Rev01