

Fluorescent Imaging Agent

Caution: For Laboratory Use. A product for research purposes only.

AngioSense® 680 EX

Product Number: NEV10054EX

DESCRIPTION: *AngioSense® 680 EX* is a fluorescent *in vivo* blood pool imaging agent. *AngioSense® 680 EX* is a near-infrared labeled fluorescent macromolecule that remains localized in the vasculature for extended periods of time and enables imaging of blood vessels and angiogenesis.

MATERIAL: (*Needs to be reconstituted*)

CONTENTS: Each vial contains 24 nmol of *AngioSense® 680 EX* as a lyophilized solid. The solution of *AngioSense® 680 EX* 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 *AngioSense® 680 EX* per mouse.

PROPERTIES: The physical properties of *AngioSense® 680 EX* can be found in **Table 1** and **Figure 1**.

STORAGE & HANDLING:

- Upon receipt, *AngioSense® 680 EX* should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *AngioSense® 680 EX* is stable for up to six months from the date of shipment.
- Once reconstituted, the PBS solution is stable up to 10 days when stored at 2-8 °C and protected from light.

IN VIVO IMAGING & APPLICATIONS:

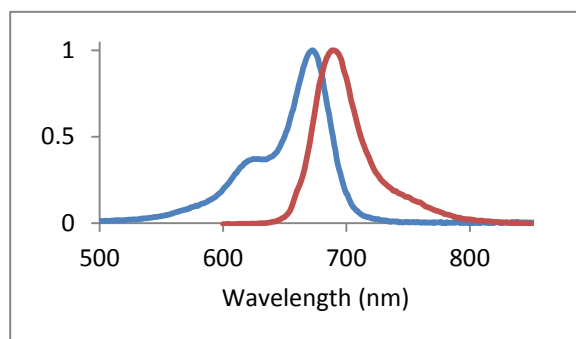
- The recommended procedure for *in vivo* imaging with *AngioSense® 680 EX* is administration via tail vein injection and imaging **24 hours post injection**.
- **Imaging in Oncology:** *AngioSense® 680 EX* can be used to study angiogenesis, as a marker for blood vessel density, in animal tumor models.
- **Imaging in Arthritis:** *AngioSense® 680 EX* can be used to characterize vascular changes and therapeutic responses associated with animal models of arthritis.

Table 1. *AngioSense® 680 EX* Properties

Property	Specification
MW	~ 70,000 g mol ⁻¹
Fluorescence ¹	
• Excitation	670 nm
• Emission	690 nm
Absorbance ¹	673 ±10 nm
Purity ²	>90%
Appearance	Blue solid

1. Absorbance, excitation, and fluorescence maxima of *AngioSense® 680 EX* in 1xPBS.
2. As determined by SE-HPLC and measuring absorbance at 680 nm.

Fig 1. Abs/Em Spectra of *AngioSense® 680 EX* in 1xPBS



NOTES:

- *PerkinElmer's AngioSense® 680 EX* is intended for research purposes only and is not for human use. It must be used by or directly under the supervision of a technically qualified individual experienced in handling potentially hazardous materials. Please read the Material Safety Data Sheet (MSDS) provided for this product.
- Several of *PerkinElmer* products and product applications are covered by U.S. and foreign patents and patents pending. Our products are not available for resale or other commercial uses without a specific agreement with *PerkinElmer*.

All other trademarks are the property of their respective owners.