

Fluorescent Imaging Agent

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

TLectinSense™ 680 Product Number: NEV10060

DESCRIPTION: *TLectinSense™ 680* is a fluorescent *in vivo* endothelial cell imaging agent. *TLectinSense™ 680* is a near-infrared labeled fluorescent macromolecule that targets the vasculature and enables imaging of blood vessels and angiogenesis.

MATERIAL: (*Needs to be reconstituted*)

CONTENTS: Each vial contains 24 nmol of *TLectinSense™ 680* as a lyophilized solid. The 1XPBS solution of *TLectinSense™ 680* has been filtered through a 0.2 µm filter prior to lyophilization. Upon reconstitution with 1.2 mL of DISTILLED WATER, this material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/100 µL of *TLectinSense™ 680* per mouse. The lyophilized formulation contains salt. When reconstituted with the recommended amount of water the salt concentration will be equivalent to 1XPBS.

PROPERTIES: The physical properties of *TLectinSense™ 680* can be found in **Table 1** and **Figure 1**

STORAGE & HANDLING:

- Upon receipt, *TLectinSense™ 680* should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.**
- When stored and handled properly, *TLectinSense™ 680* is stable for 3 months from the date of shipment.
- Once reconstituted, the solution is stable up to 10 days when stored at 2-8 °C and protected from light.

IN VIVO IMAGING AND APPLICATIONS:

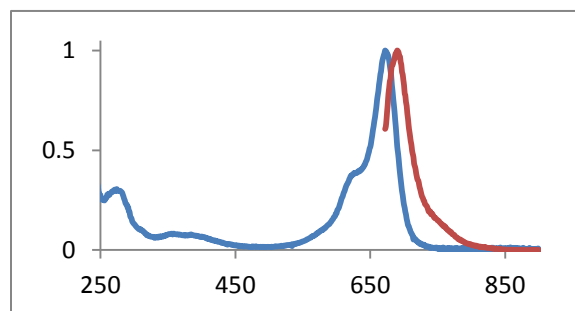
- The recommended procedure for *in vivo* imaging with *TLectinSense™ 680* is intravenous administration and imaging **6 hours post injection.**
- **Imaging in matrigel plugs and tumor models:** *TLectinSense™ 680* can be used to study angiogenesis and blood vessel density, in matrigel plugs and animal tumor models.

Table 1. *TLectinSense™ 680* Properties

Property	Specification
MW	~ 72,000 g mol ⁻¹
Fluorescence ¹	
• Excitation	670 nm
• Emission	690 nm
Absorbance ¹	675 ± 10nm
Purity ²	>95%
Appearance	Blue solid

1. Absorbance, excitation, and fluorescence maxima in PBS.
2. As determined by SE-HPLC and measuring absorbance at 675 nm.

Fig 1. Abs/Em Spectra of *TLectinSense™ 680* in 1xPBS



NOTES:

- *PerkinElmer's TLectinSense™ 680* 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*.

PerkinElmer, Inc.
549 Albany Street
Boston, MA 02118 USA
P: (800) 762-4000 or (+1) 203-925-4602
www.perkinelmer.com/enradiochemicals

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc.
All other trademarks are the property of their respective owners.

