

# Fluorescent Imaging Agent

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

## BacteriSense™ 645

Product Number: NEV10080

### DESCRIPTION

*BacteriSense™ 645* is a near-infrared labeled fluorescent *in vivo* imaging agent that targets Gram positive and Gram negative bacteria.

### CONTENTS

- Each vial contains 60 nmol of *BacteriSense™ 645* as a lyophilized solid.
- The aqueous solution of *BacteriSense™ 645* 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 5 nmol/100 µL of *BacteriSense™ 645* per mouse.

### STORAGE & HANDLING

- Upon receipt, *BacteriSense™ 645* should be IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT.
- When stored and handled properly, *BacteriSense™ 645* is stable for up to three months from the date of shipment.
- To prevent material losses, vials should be centrifuged before opening. Measures should be taken to reduce dispersion due to static charge. Use of a static gun is recommended.
- Once reconstituted, the solution is stable up to 10 days when stored at 2-8 °C.

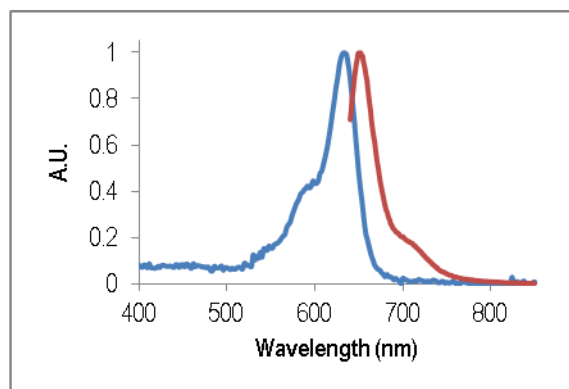
### IN VIVO IMAGING AND APPLICATIONS

- The recommended procedure for *in vivo* imaging with *BacteriSense™ 645* is intravenous administration and imaging 1 hour post injection.
- *BacteriSense™ 645* can be used to study superficial and deep tissue bacterial infections due to a variety of bacteria including *Escherichia coli* and *Staphylococcus epidermidis*, as well as the effect of antibiotic treatments.

### PHYSICAL AND SPECTRAL PROPERTIES

Property	Specification
MW	1,010 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	635 nm
• Emission	656 nm
Absorbance <sup>1</sup>	635 ±10nm
Purity <sup>2</sup>	>95%
Appearance	Blue solid

1. Absorbance and fluorescence maxima in 1 x PBS.
2. As determined by SE-HPLC; measuring absorbance at 635 nm.



Absorbance and fluorescence emission spectra in 1x PBS.

## NOTES

- *PerkinElmer's BacteriSense™ 645* 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/nenradiochemicals](http://www.perkinelmer.com/nenradiochemicals)

**For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://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.

