

# Fluorescent Imaging Agent

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

## Genhance™ 750

**Product Number: NEV10118**

**DESCRIPTION:** *Genhance*™ 750 is the acid form of VivoTag® 680 that can be used as a control or a fluorescent *in vivo* blood pool imaging agent.

Upon dilution, a 1 mg package of *Genhance* 750 provides sufficient reagents for imaging approximately 20 mice (weighing ~25 grams each) when using the recommended dose of 0.05mg/100 µL 1xPBS.

### MATERIAL

**CONTENTS:** *Genhance* 750: Available as a solid in 1 mg vials.

**PROPERTIES:** The physical characteristics of *Genhance* 750 can be found in **Table 1** and **Figure 1**.

### STORAGE & HANDLING:

- Upon receipt, *Genhance* 750 should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *Genhance* 750 is stable for up to twenty four months.
- Allow *Genhance* 750 imaging agent to equilibrate to room temperature before opening.

### IN VIVO IMAGING AND APPLICATIONS:

- Dilute *Genhance* 750 to a concentration of 0.5mg/mL with sterile 1X PBS.
- The generally recommended procedure for *in vivo* imaging with *Genhance* 750 is administration via tail vein injection and imaging **30 minutes post injection**.
- *Genhance* 750 can be used as a control during imaging studies and as a blood pool imaging agent.

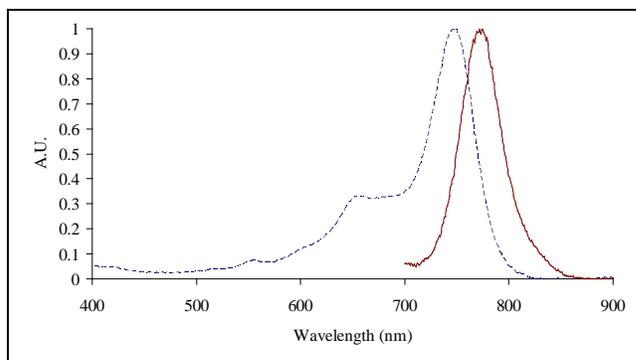
**Table 1. *Genhance* 750 Characteristics**

| Property                  | Specification                            |
|---------------------------|--|
| MW                        | 1086 g mol <sup>-1</sup>                 |
| Fluorescence <sup>1</sup> |  |
| • Excitation              | 750 ±5 nm                                |
| • Emission                | 775 ±5 nm                                |
| Absorbance <sup>1</sup>   | 750 ±5nm                                 |
| Extinction                | ~240,000M <sup>-1</sup> cm <sup>-1</sup> |
| Purity <sup>2</sup>       | >90%                                     |
| Appearance                | Blue Solid                               |

1. Absorbance and fluorescence maxima determined in 1x PBS.

2. As determined by RP-HPLC and measuring absorbance at 750 nm.

\* MW as Bis TEA Salt



**Figure 1.** Absorbance and fluorescence emission spectra of

## NOTES:

- *PerkinElmer's Genhance 750* 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's* 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 from *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.

