

Fluorescent Imaging Agent

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

Genhance™ 750

Product Number: NEV10177

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

Upon dilution, a 1mg 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: Solid 5 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 ⁻¹
Fluorescence ¹	
• Excitation	750 ± 5 nm
• Emission	775 ± 5 nm
Absorbance ¹	750 ± 5 nm
Extinction	~240,000 M ⁻¹ cm ⁻¹
Purity ²	>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

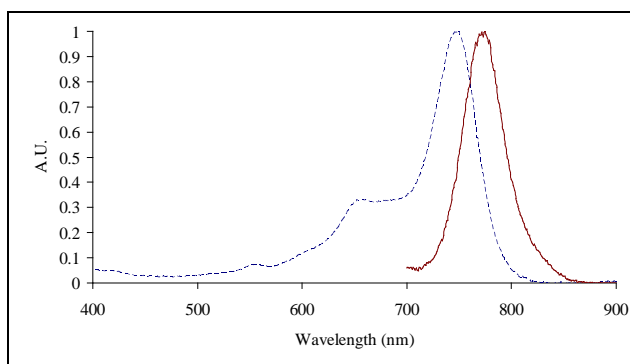


Figure 1. Absorbance and fluorescence emission spectra of

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 with *PerkinElmer*.

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

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.

