

# Fluorescent Imaging Agent

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

## PSA 750 FAST™

Product Number: NEV11125

### DESCRIPTION

PSA 750 FAST™ is a prostate-specific antigen (PSA) activatable agent that is optically silent upon injection and produces fluorescent signal after cleavage by enzymatically active PSA. Activation only occurs in the presence of active PSA. PSA 750 FAST™ may be used to monitor the progression of prostate tumors or to evaluate the potential therapeutic efficacy of drugs targeting the underlying mechanisms involved in these diseases.

### CONTENTS

- Each vial contains 24 nmol of PSA 750 FAST™ in dry solid form. PSA 750 FAST™ has been filtered through a 0.2µm filter prior to drying.
- Reconstitute PSA 750 FAST™ with 1.2 mL of 1X PBS before administering to animals.
- The packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol (in 100 µL of PBS) of PSA 750 FAST™ per mouse.

### STORAGE & HANDLING

- Upon receipt, PSA 750 FAST™ should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, PSA 750 FAST™ in its dry solid form is stable for up to six months.
- Before opening the vial check to ensure that all of the solid material is at the bottom of the vial.
- After reconstituting with 1X PBS, gently swirl the solution to ensure that the solid is fully in solution.
- **Once reconstituted, PSA 750 FAST™ is stable for up to 7 days when stored at 2-8 °C and protected from light.**
- Allow PSA 750 FAST™ imaging agent to equilibrate to room temperature before introducing into animals.

### IN VIVO IMAGING AND APPLICATIONS

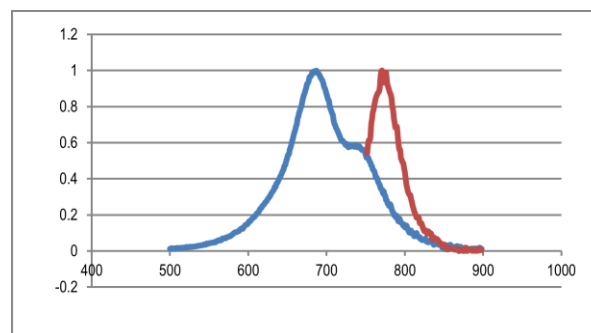
- The generally recommended procedure for *in vivo* imaging with PSA 750 FAST™ is administration *via* intravenous injection and imaging **6-24 hours post injection**.
- PSA 750 FAST™ will clear from tissues after approximately 50 hours. Repeat injection and imaging may be performed every 7 days for longitudinal studies. It is recommended that a pre-injection baseline image be taken prior to reinjection and imaging.
- PSA 750 FAST™ enables imaging of PSA positive tumors such as LNCaP implanted in male nude mice.

### PHYSICAL AND SPECTRAL PROPERTIES

Property	Specification
MW	~42,500 g mol <sup>-1</sup>
Fluorescence emission <sup>1</sup>	770 nm
Absorbance <sup>1</sup>	687 nm (quenched) 750 nm (activated)
Purity <sup>2</sup>	>95 %
Appearance	Blue solid

1. Absorbance and fluorescence maxima in 1x PBS.

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



Normalized absorbance (blue) and fluorescence emission (red) spectra of PSA 750 FAST™ in 1x PBS

## NOTES

- *PSA 750 FAST™* 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.

