

Fluorescent Imaging Agent

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

ProSense® 750 EX

Product Number: NEV10001EX

DESCRIPTION: *ProSense*®750 EX is a protease activatable fluorescent *in vivo* imaging agent that is activated by key disease associated proteases such as Cathepsin B, L, S and Plasmin. *ProSense*®750 EX is optically silent in its unactivated state and becomes highly fluorescent following protease-mediated activation.

MATERIAL: (*Needs to be reconstituted*)

CONTENTS: Each vial contains 24 nmol (fluorophore) of *ProSense*®750 EX as a lyophilized solid. The solution of *ProSense*®750 EX 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 2 nmol/100 µL 1xPBS of *ProSense*®750 EX per mouse.

PROPERTIES: The physical properties of *ProSense*®750 EX can be found in **Table 1** and **Figure 1**.

STORAGE & HANDLING:

- Upon receipt, *ProSense*®750 EX should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *ProSense*®750 EX is stable for up to six months from the date of shipment.
- Once reconstituted, the PBS solution is stable up to 10 days when stored at 2-8 °C.

IN VIVO IMAGING & APPLICATIONS:

- The recommended procedure for *in vivo* imaging with *ProSense*®750 EX is administration via tail vein injection and imaging **24 hours post injection**.
- **Imaging in Arthritis:** *ProSense*®750 EX can be used as a marker for disease progression and therapeutic response in animal models of arthritis.
- **Imaging in Oncology:** *ProSense*®750 EX can be used as a marker for disease progression in animal tumor models.

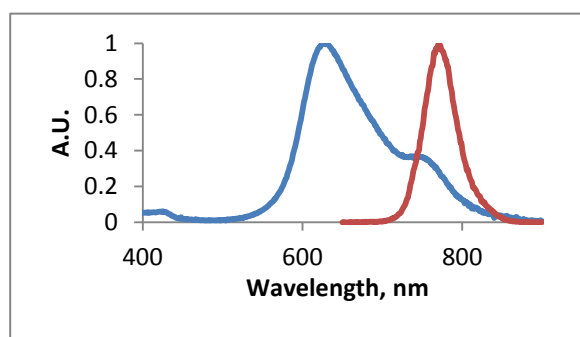
Table 1. *ProSense*®750 EX Properties

Property	Specification
MW	~ 450,000 g mol ⁻¹
Fluorescence ¹	
• Excitation	750 nm
• Emission	770 nm
Absorbance ¹	630 ±10nm
Purity ²	>95%
Appearance	Blue solid

1. Absorbance, excitation, and fluorescence maxima of *ProSense*®750 EX in 1xPBS.

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

Fig 1. Abs/Em Spectra of *ProSense*®750 EX in 1xPBS



NOTES:

- *PerkinElmer's ProSense[®] 750 EX* 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

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.

