

# Fluorescent Imaging Agent

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

## ProSense® 680

Product Number: NEV10003

**DESCRIPTION:** *ProSense*® 680 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* 680 is optically silent in its unactivated state and becomes highly fluorescent following protease-mediated activation.

**MATERIAL:** (*Needs to be reconstituted*)

**CONTENTS:** Each vial contains 20 nmol of *ProSense* 680 at a concentration of 20 nmol/150 µL, in 1xPBS. The *ProSense* 680 solution has been filtered through a 0.2 µm filter. Upon dilution with 1350 ul of 1 x PBS, this material provides sufficient reagents for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/150 µL 1xPBS of *ProSense* 680 per mouse.

**PROPERTIES:** The physical properties of *ProSense* 680 can be found in **Table 1 and Figure 1.**

### STORAGE & HANDLING:

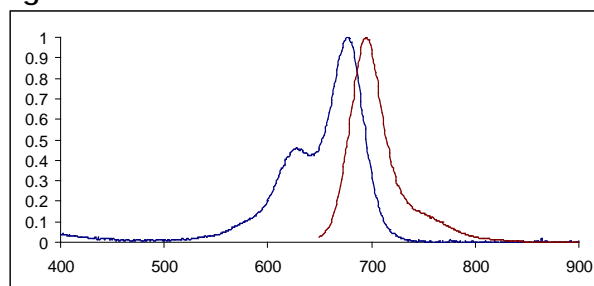
- Upon receipt, *ProSense* 680 should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.**
- When stored and handled properly, *ProSense* 680 is stable for up to twelve months.
- Allow *ProSense* 680 imaging agent to equilibrate to room temperature before injecting into animals.

### IN VIVO IMAGING & APPLICATIONS:

- The recommended procedure for *in vivo* imaging with *ProSense* 680 is administration via tail vein injection and imaging **24 hours post tail vein injection.**
- **Imaging in Arthritis:** *ProSense* 680 can be used as a marker for disease progression and therapeutic response in animal models of arthritis.
- **Imaging in Oncology:** *ProSense* 680 can be used as a marker for disease progression in animal tumor model

Property	Specification
MW	~400,000 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	<ul style="list-style-type: none"> <li>• Excitation 680 ±10 nm</li> <li>• Emission 700 ±10 nm</li> </ul>
Absorbance <sup>1</sup>	680 ±10 nm
Purity <sup>2</sup>	>95%
Appearance	Clear blue solution

Fig 1.



Absorbance and fluorescence emission spectra in 1x PBS.

## SELECTED REFERENCES:

- Weissleder, R., Tung, C.H., Mahmood, U., Bogdanov, A. In vivo imaging of tumors with protease-activated near-infrared fluorescent probes. *Nature Biotechnology* **17**, 375-378 (1999)

## NOTES:

- *PerkinElmer's ProSense 680* 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](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.

