

# Fluorescent Imaging Agent

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

## MMPSense™ 680

**Product Number: NEV10126**

**DESCRIPTION:** *MMPsense™ 680* is a protease activatable fluorescent *in vivo* imaging agent that is activated by key matrix metalloproteinases (MMPs) including MMP-2, -3, -9 and -13. *MMPsense 680* is optically silent in its unactivated state and becomes highly fluorescent following protease-mediated activation.

**MATERIAL:** (Ready to use: No dilution required)

**CONTENTS:** Each vial contains 20 nmol of *MMPsense 680* at a concentration of 20 nmol/1.5 mL in 1xPBS. The *MMPsense 680* solution has been filtered through a 0.2 µm filter. This package provides sufficient reagents for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol/150 µL 1xPBS of *MMPsense 680* per mouse.

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

### STORAGE & HANDLING:

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

### IN VIVO IMAGING AND APPLICATIONS:

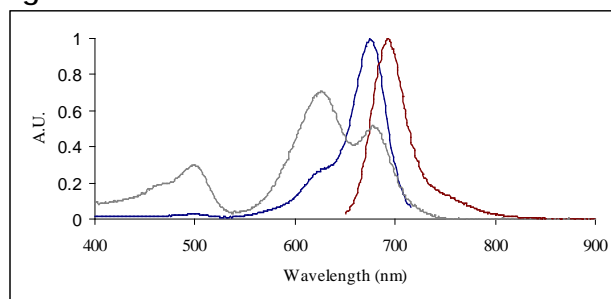
- The recommended procedure for *in vivo* imaging with *MMPsense 680* is administration via tail vein injection and imaging **24 hours post tail vein injection**.
- **Imaging in Oncology:** *MMPsense 680* can be used as a marker for disease progression in animal tumor models.

**Table 1.** *MMPsense 680* Properties

Property	Specification
MW	~ 450,000 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	680 ± 10 nm
• Emission	700 ± 10nm
Absorbance <sup>1</sup>	680 ± 10nm
Purity <sup>2</sup>	>95%
Appearance	Clear Blue Solution

1. Absorbance and fluorescence maxima of *MMPsense 680* in PBS.
2. As determined by SE-HPLC and measuring absorbance at 680 nm.

**Fig 1.**



## NOTES:

- *PerkinElmer's MMPsense 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.
- 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.

