

# Fluorescent Imaging Agent

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

## FolateRSense™ 680

Product Number: NEV10040

**DESCRIPTION:** *FolateRSense™ 680* is a folate receptor targeted fluorescent *in vivo* imaging agent that can be used to image folate receptor expression in tumors.

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

**CONTENTS:** Each vial contains 24 nmol of *FolateRSense™ 680* in dry solid form. Reconstitute *FolateRSense™ 680* with 1.2 mL of 1 x PBS before injecting into animals. Each vial of packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol (100 µL) of *FolateRSense™ 680* per mouse.

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

### STORAGE & HANDLING:

- Upon receipt, *FolateRSense™ 680* should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *FolateRSense™ 680* is stable for 3 months from the date of shipment.
- Before opening the vial check to ensure that all of the solid material is at the bottom of the vial.
- After reconstituting with PBS, gently swirl the solution to ensure that the solid is fully in solution.
- **Once reconstituted with 1 x PBS, the solution is stable up to 10 days when stored at 2-8 °C and protected from light.**

### IN VIVO IMAGING AND APPLICATIONS:

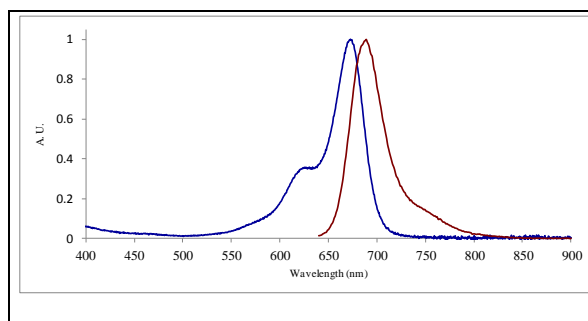
- The generally recommended procedure for *in vivo* imaging with *FolateRSense™ 680* is administration via intravenous injection and imaging **6-24 hours post injection**.

**Table 1.** *FolateRSense™ 680* Properties

Property	Specification
MW	1605.7 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	
• Excitation	670 nm
• Emission	690 nm
Absorbance <sup>1</sup>	670 nm
Purity <sup>2</sup>	>90%
Appearance	Blue solid

1. Absorbance and fluorescence maxima in 1x PBS.  
2. Determined by RP-HPLC and measuring absorbance at 670 nm.

**Fig 1.** Abs/Em Spectra of *FolateRSense™ 680* in 1xPBS



Normalized absorbance (blue) and fluorescence emission (red) spectra in 1x PBS.

## NOTES:

- *FolateRSense*<sup>TM</sup> 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*.