

Fluorescent Imaging Agent

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

Superhance® 680

Product Number: NEV10116

DESCRIPTION: *Superhance*TM 680 is a small molecule fluorescent *in vivo* blood pool imaging agent. *Superhance* 680 enables imaging of circulation, blood vessels, vasculature, vascular leak, including that associated with early oncologic and ophthalmologic lesions.

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

CONTENTS: Each vial contains 4 nmoles of *Superhance* 680 in 1500 ul at a concentration of 4 nmol/150 µL, in 1xPBS. The *Superhance* 680 solution has been filtered through a 0.2 µm filter. This material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 4 nmoles of *Superhance* 680 per mouse.

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

STORAGE & HANDLING:

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

IN VIVO IMAGING AND APPLICATIONS:

- The generally recommended procedure for *in vivo* imaging with *Superhance* 680 is administration via tail vein injection and imaging **0.5 – 24 hours post injection**.
- *Superhance* 680 enables imaging of circulation, blood vessels, vasculature, vascular leak, including that associated with early oncologic and ophthalmologic lesions.

Table 1. Superhance 680 Characteristics

Property	Specification
MW	~ 1540 g mol ⁻¹
Fluorescence ¹	
• Excitation	675 ± 5 nm
• Emission	692 ± 5 nm
Absorbance ¹	675 ± 5 nm
Purity ²	>95%
Appearance	Clear blue solution

1. Absorbance and fluorescence maxima of *Superhance* 680 in 1x PBS.

2. As determined by RP-HPLC and measuring absorbance at 675 nm.

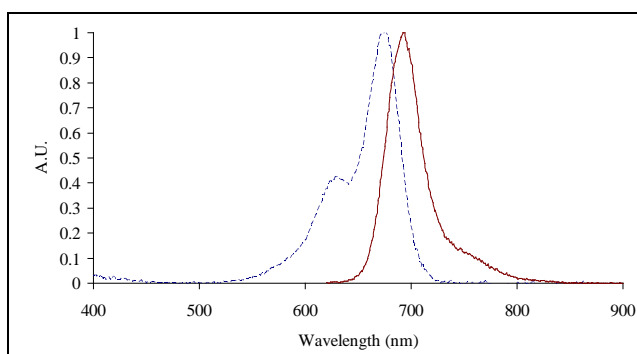


Figure 1. Absorbance and fluorescence emission spectra of *Superhance* 680 in 1x PBS.

NOTES:

- *PerkinElmer's Superhance 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 from *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.

