

Fluorescent Imaging Agent

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

IntegriSense™ 680

Product Number: NEV10645

DESCRIPTION: *IntegriSense™ 680* is a targeted fluorescence imaging agent comprising a potent, selective non-peptide small molecule integrin $\alpha_v\beta_3$ antagonist and an NIR fluorochrome. This agent has been developed to enable *in vivo* visualization and quantification of integrin $\alpha_v\beta_3$ expression in neovasculature as well as in tumor cells, to monitor tumor angiogenesis, growth and treatment efficacy. Half-life in tissue of IntegriSense 680 signal is >21 days.

MATERIAL (Needs to be reconstituted)

CONTENTS: Each vial contains 24 nmol of *IntegriSense 680* lyophilized solid. The *IntegriSense 680* solution has been filtered through a 0.2 μm filter prior to lyophilization. Reconstitute *IntegriSense 680* with 1.2 mL of 1 x PBS before injecting into animals. The 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 *IntegriSense 680* per mouse.

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

STORAGE & HANDLING:

- Upon receipt, *IntegriSense 680* should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT.**
- When stored and handled properly, *IntegriSense 680* is stable for up to twelve months in the lyophilized form.
- Before opening the vial check to ensure that the lyophilized powder is present at the bottom of the vial.
- After reconstituting with PBS, gently swirl the solution to ensure that the lyophilized powder 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.**

Table 1. *IntegriSense 680* Characteristics

Property	Specification
MW	1432 g mol ⁻¹
Fluorescence ¹	
• Excitation	675 nm
• Emission	693 nm
Absorbance	675 \pm 5 nm
Purity ²	>95 %
Appearance	Dark blue solid

1. Absorbance and fluorescence maxima of *IntegriSense 680* in 1x PBS.
2. As determined by RP-HPLC and measuring absorbance at 675 nm.

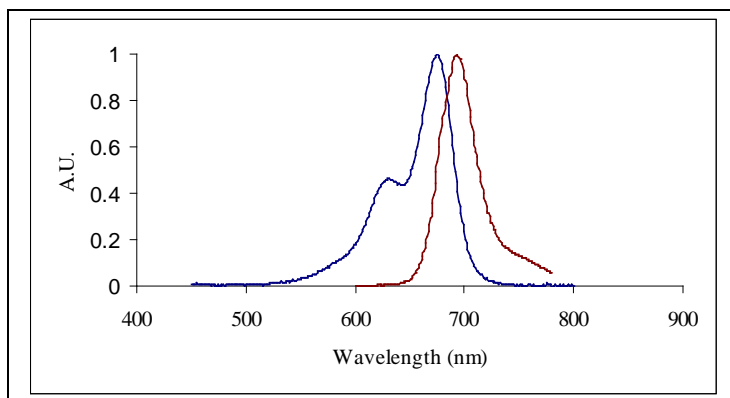


Figure 1. Absorbance and fluorescence emission spectra of *IntegriSense 680* in 1x PBS.

IN VIVO IMAGING AND APPLICATIONS:

- The generally recommended procedure for *in vivo* imaging with *IntegriSense 680* is administration via intravenous injection and imaging **3 – 48 hours post injection**.
- Imaging at earlier time points (~3 hrs) is recommended when imaging the vasculature
- Imaging at later time points (24-48 hrs) is recommended when imaging tumors to reduce background.
- *IntegriSense 680* enables imaging of neovasculature and tumors in a range of oncology applications.

NOTES:

- *PerkinElmer's IntegriSense 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.

