

Fluorescent Imaging Agent

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

AngioSPARK[®] 750

Product Number: NEV10150

DESCRIPTION: NanoSPARKS[™] are a new family of highly fluorescent near infrared nanoparticles specifically designed for *in vivo* imaging. NanoSPARKS contain an iron oxide core that is coated to specifically produce a functionalized biocompatible product. AngioSPARK[®] 750 comprises pegylated fluorescent nanoparticles that remain localized in the vasculature for extended periods of time and enable imaging of blood vessels and angiogenesis.

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

CONTENTS: Each vial contains 500 µL of AngioSPARK 750 in 1xPBS. The AngioSPARK 750 solution has been filtered through a 0.2 µm filter.

This material provides sufficient reagent for imaging approximately 5 mice (weighing ~25 grams each) when using the recommended injection volume of 100 µL of AngioSPARK 750 per mouse.

PROPERTIES: The physical characteristics of AngioSPARK 750 can be found in **Table 1** and **Figure 1**.

STORAGE & HANDLING:

- Upon receipt, AngioSPARK 750 should be **IMMEDIATELY STORED AT 2-8 °C AND PROTECTED FROM LIGHT. DO NOT FREEZE.**
- When stored and handled properly, AngioSPARK 750 is stable for up to 12 months.
- Allow AngioSPARK 750 to equilibrate to room temperature before use.

IN VIVO IMAGING APPLICATION:

- The recommended procedure for analysis by intravital microscopy imaging using AngioSPARK 750 is administration via tail vein injection and imaging **0 – 4 hour post injection.**
- AngioSPARK 750 can also be used at later time points (>12 hours) to image vascular leak in tissue site of inflammation and cancer. AngioSPARK 750 can be imaged within the interstitium for up to 24 hours post tail vein injection.

NOTES:

Table 1. AngioSPARK 750 Characteristics

Property	Specification
Particle Size	20-50nm
Fluorescence ¹	
• Excitation	750 ± 5 nm
• Emission	775 ± 5 nm
Extinction ¹	4 X 10 ⁶ ± 2 X 10 ⁶ M ⁻¹ cm ⁻¹
Appearance	Greenish-brown solution

1. Absorbance and fluorescence maxima of AngioSPARK 750 in 1x PBS

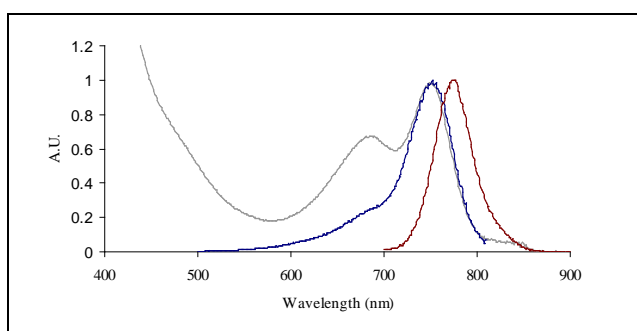


Figure 1. Absorbance, excitation and fluorescence emission spectra of AngioSPARK 750 in 1x PBS.

- *PerkinElmer's* NanoSPARK's products are intended for research purposes only and are 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.

