

# NIR Fluorochrome Label

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

## VivoTag® 680 XL *IN VIVO* NIR FLUOROCHROME LABEL

Product Number: NEV11119

**DESCRIPTION:** VivoTag® 680 XL is an amine reactive near-infrared (NIR) fluorochrome (NHS ester) for labeling biomolecules for *in vivo* imaging applications.

### MATERIAL

**CONTENTS:** *VivoTag 680 XL*: Available in 1 mg or 5 mg vials.

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

### STORAGE & HANDLING:

- Upon receipt, VivoTag 680 XL should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *VivoTag 680 XL* is stable for up to three months.
- Allow *VivoTag 680 XL* to equilibrate to room temperature before using.

### IN VIVO IMAGING AND APPLICATIONS:

- Optimized *in vivo* NIR label for labeling target molecules including peptides, small molecules, macromolecules and nanoparticles.
- Highly hydrophilic and suitable for multiple dye loading.

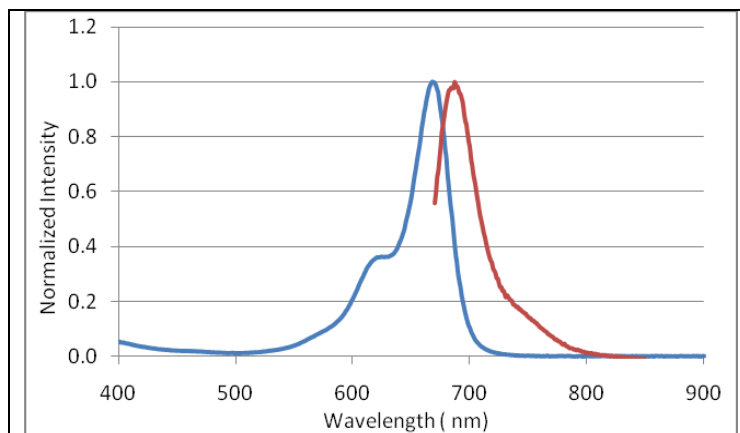
**Table 1.** *VivoTag 680 XL* Characteristics

Property	Specification
MW	1856 g mol <sup>-1</sup>
Fluorescence <sup>1</sup>	<ul style="list-style-type: none"> <li>• Excitation 665 ± 5 nm</li> <li>• Emission 688 ± 5 nm</li> </ul>
Absorbance <sup>1</sup>	668 ± 3 nm
Extinction	~210,000 M <sup>-1</sup> cm <sup>-1</sup>
Purity <sup>2</sup>	>95%
Appearance	Blue solid

1. Absorbance and fluorescence maxima determined in 1x PBS.

2. Reactive ester as determined by RP-HPLC and measuring absorbance at 670 nm.

\*MW as Penta TEA Salt



Absorbance and fluorescence emission spectra of *VivoTag 680 XL* in 1x PBS.

## NOTES:

- *PerkinElmer's VivoTag 680 XL* 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](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.

