

NIR Fluorochrome Label

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

VivoTag-S[®] 680

Product Number: NEV10122

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

MATERIAL

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

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

STORAGE & HANDLING:

- Upon receipt, *VivoTag-S 680* should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *VivoTag-S 680* is stable for up to twenty four months.
- Allow *VivoTag-S 680* 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.
- Labeling of target molecules that have multiple amine reactive sites with > 1 fluorochrome results in a blue shift in the absorbance spectra as shown in **Fig. 1**.

Table 1. *VivoTag-S 680* Characteristics

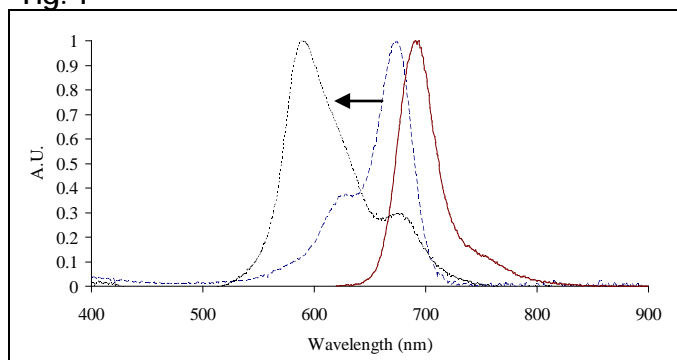
Property	Specification
MW	1439 g mol ⁻¹
Fluorescence ¹	<ul style="list-style-type: none"> • Excitation 673 ± 5 nm • Emission 691 ± 5 nm
Absorbance ¹	673 ± 5 nm
Extinction	~230,000 M ⁻¹ cm ⁻¹
Purity ²	>90 %
Appearance	Blue Solid

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

2. Reactive ester as determined by RP-HPLC and measuring absorbance at 675 nm.

* MW as Tris TEA salt

Fig. 1



Absorbance and fluorescence emission spectra of *VivoTag-S 680* in 1x PBS. Arrow indicates shift in absorbance of conjugate with multiple fluorochromes.

NOTES:

- *PerkinElmer's VivoTag-S 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.

