

NIR Fluorochrome Label

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

VivoTag® 800 *IN VIVO* NIR FLUOROCHROME LABEL

Product Number: NEV11107

DESCRIPTION: *VivoTag*® 800 is an amine reactive (NHS Ester) near-infrared (NIR) fluorochrome for labeling biomolecules for *in vivo* imaging applications.

MATERIAL

CONTENTS: *VivoTag* 800: Available in 1 mg or 5 mg vials.

PROPERTIES: The physical characteristics of *VivoTag* 800 can be found in **Table 1** and **Figure 1**.

STORAGE & HANDLING:

- Upon receipt, *VivoTag* 800 should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *VivoTag* 800 is stable for up to three months.
- Allow *VivoTag* 800 to equilibrate to room temperature before using.
- **Recommended solvents for reconstitution:**
Aqueous buffers pH 6.0 -7.0
Organic solvents: DMSO, DMF
- **Recommended solvents for performing conjugation reactions:** Aqueous buffers in the pH 7.5 - 8.5 range (borate or phosphate buffer) or DMSO or DMF
- Carbonate buffers are not recommended for conjugation reactions

IN VIVO IMAGING AND APPLICATIONS:

- Optimized *in vivo* NIR label for labeling target molecules including peptides, small molecules, macromolecules and nanoparticles.

Table 1. *VivoTag* 800 Characteristics

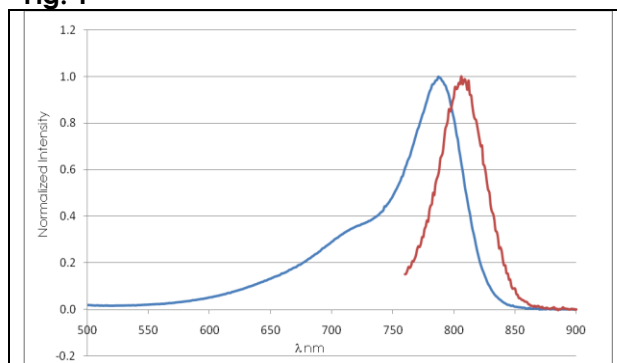
Property	Specification
MW	1464 g mol ⁻¹
Fluorescence ¹	<ul style="list-style-type: none"> • Excitation 785 nm • Emission 810 nm
Absorbance ¹	787 ± 5 nm
Extinction	~200,000 M ⁻¹ cm ⁻¹
Purity ²	>90%
Appearance	Green solid

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

2. Reactive ester as determined by RP-HPLC and measuring absorbance at 790 nm.

* MW as tris triethylamine salt

Fig. 1



Absorbance and fluorescence emission spectra of *VivoTag* 800 in 1x PBS

NOTES:

- *PerkinElmer's VivoTag 800* 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.

