

Fluorescent Imaging Agent

Caution: For laboratory use only. This product is intended for animal research only and not for use in humans.

Name: Transferrin-Vivo 750

Product Number: NEV10091

DESCRIPTION

- ▶ *Transferrin-Vivo 750* is a fluorescently-labeled protein designed to bind to transferrin receptors over-expressed in cancer cell with rapid proliferation.

CONTENTS

- Each vial contains 24 nmol of *Transferrin-Vivo 750* in dry solid form. *Transferrin-Vivo 750* has been filtered through a 0.2µm filter prior to drying.
- Reconstitute *Transferrin-Vivo 750* with 1.2 mL of 1X PBS before administering to animals.
- The packaged material provides sufficient reagent for imaging approximately 10 mice (weighing ~25 grams each) when using the recommended dose of 2 nmol (in 100 µL of PBS) of *Transferrin-Vivo 750* per mouse.

STORAGE & HANDLING

- Upon receipt, *Transferrin-Vivo 750* should be **IMMEDIATELY STORED AT -20 °C AND PROTECTED FROM LIGHT**.
- When stored and handled properly, *Transferrin-Vivo 750* in its dry solid form is stable for up to three months.
- Before opening the vial check to ensure that all of the solid material is at the bottom of the vial.
- After reconstituting with 1X PBS, gently swirl the solution to ensure that the solid is fully in solution.
- **Once reconstituted, *Transferrin-Vivo 750* is stable for up to 14 days when stored at 2-8 °C and protected from light.**
- Allow *Transferrin-Vivo 750* imaging agent to equilibrate to room temperature before introducing into animals.

IN VIVO IMAGING AND APPLICATIONS

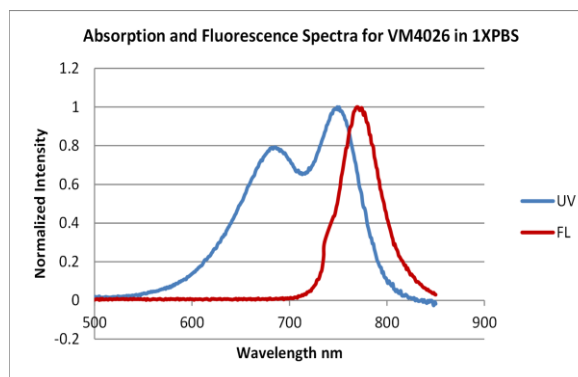
- The generally recommended procedure for *in vivo* imaging with *Transferrin-Vivo 750* is administration via intravenous injection and imaging 24 hours post injection.
- The tissue half-life of *Transferrin-Vivo 750* is approximately 24 hours. Repeat injection and imaging may be performed every 4 days for longitudinal studies. It is recommended that a pre-injection baseline image be taken prior to re-injection and imaging if repeat injections are to be performed less than 1 week apart.
- *Transferrin-Vivo 750* enables imaging of HT-29 tumors implanted subcutaneously in nude mice.

PHYSICAL AND SPECTRAL PROPERTIES

Property	Specification
MW	~106,000 g mol ⁻¹
Fluorescence emission ¹	770 nm
Absorbance ¹	750 nm
Purity ²	>90 %
Appearance	Blue solid

1. Absorbance and fluorescence maxima in 1x PBS.

2. As determined by RP-HPLC; measuring absorbance at 750 nm.



Normalized absorbance (blue) and fluorescence emission (red) spectra of *Transferrin-Vivo 750* in 1x PBS

NOTES

- Caution: For laboratory use only. This product is intended for animal research only and not for use in humans. 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.

