

XenoLight DiR

Research Use Only. Not for use in diagnostic procedures.

XenoLight DiR

Product Number: 125964

DESCRIPTION:

XenoLight DiR is a lipophilic, near infrared fluorescent cyanine dye ideal for staining the cytoplasmic membrane. The two long 18-carbon chains insert into the cell membrane, resulting in specific and stable cell staining with negligible dye transfer between cells.

XenoLight DiR can be used for non-invasive imaging of cell homing (T cells, stem cells etc.) *in vivo*. The near infrared property of this dye makes it ideal for *in vivo* imaging because of significantly reduced auto-fluorescence from the animal at higher wavelengths.

CONTENTS: 25mg of XenoLight DiR (DiIC18(7) or 1,1'-dioctadecyltetramethyl indotricarbocyanine Iodide)

PROPERTIES: The physical properties of XenoLight DiR can be found in **Table 1**.

STORAGE & HANDLING:

- Store at $\leq 4^{\circ}\text{C}$ and protect from light.

IN VIVO IMAGING & APPLICATIONS:

- A stock solution of the dye can be made in ethanol
- Cell staining can be effected using the dye at 1-10 μM concentration and 10-20 minute incubation time.

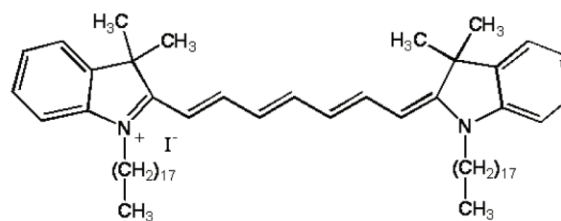
Key Applications:

Cell Staining

XenoLight DiR can be applied to fluorescence staining of primary cells, such as embryonic stem cells, bone marrow derived stem cells, adipose derived stem cells, lymphocytes and erythrocytes, which will enable fluorescence detection of stained cells and their *in vivo* distribution. Since DiR has excitation and emission maxima in the NIR range, fluorescence detection of DiR stained cells will have less interference from auto-fluorescent tissue background, resulting in high sensitivity of detection.

Table 1. XenoLight DiR Properties

Property	Specification
Molecular Weight	1013.4
Fluorescence <ul style="list-style-type: none"> • Absorption • Emission 	748 nm (in MeOH) 780 nm (in MeOH)
IVIS Filter Set	710 ex/760 em
Extinction	270,000 (in MeOH)
Color and Form	Blue-green oily solid
Solubility	Soluble in ethanol, DMF or DMSO



In Vivo Imaging of DiR Stained Spleen T-cell Distribution

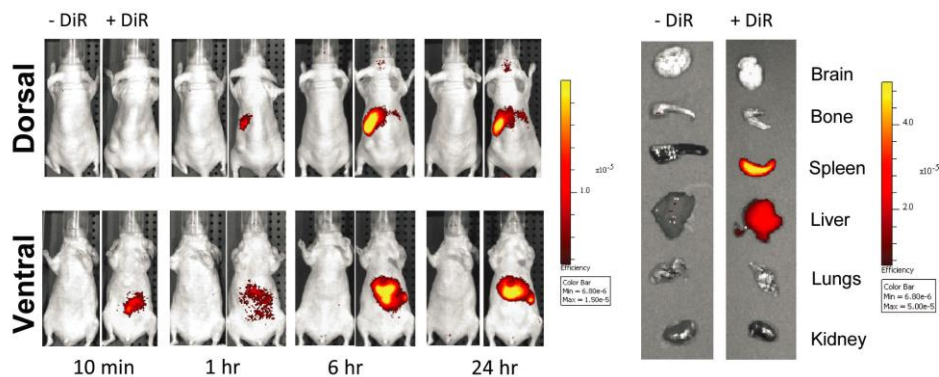


Figure 1. XenoLight DiR stock was prepared by dissolving 25 mg in 3 mL ethanol. Working solution of 320 $\mu\text{g}/\text{mL}$ was prepared by diluting 199 μL of stock solution in 5 mL PBS. T-cells isolated from the spleen were incubated with 320 $\mu\text{g}/\text{mL}$ DiR. After 30 min incubation, cells were spun down for 3 min at 1000 rpm at 4 $^{\circ}\text{C}$ resulting in a blue pellet. Cells were washed twice in PBS and injected intravenously (5×10^6 cells/mouse). Control group was injected with 5×10^6 cells/mouse in PBS. Mice were imaged with IVIS Spectrum at 10 min, 1 hr, 6 hr and 24 hrs post injection. Ideal filter set for DiR imaging is 710 nm excitation and 760 nm emission. Mice were imaged dorsally as well as ventrally at all time points. Brain, bones, spleen, liver, lungs and kidneys were harvested for ex vivo imaging 24 hrs post injection. Non-invasive in vivo imaging showed the homing process of injected T cells to the liver and spleen in real time, which was confirmed by ex vivo imaging.

PRODUCT WARRANTY:

PerkinElmer warrants that this product will be stable upon shipment from PerkinElmer for a period of twelve months, provided that it has been properly stored and protected from light during this period.

NOTES:

Reference:

Kalchenko et al., Use of lipophilic near-infrared dye in whole-body optical imaging of hematopoietic cell homing. *Journal of Biomedical Optics*, September/October 2006, Vol 11(5).

Label License:

This product is provided under an Agreement between Biotium and PerkinElmer (formerly Caliper Life Sciences). The manufacture, use, sale or import of this product is subject to one or more of pending patent applications owned by Biotium. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer, where such research does not include testing, analysis or screening services for any third party in return for compensation on a per test basis. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. For information on purchasing a license to this product for purposes other than research, contact Biotium, Inc., Business Development, 3423 Investment Blvd, Suite 8, Hayward, CA 94545, Tel: 510-265-1027, Fax: 510-265-1352.

Distributed by:
PerkinElmer, Inc.
549 Albany Street
Boston, MA 02118 USA
P: (800) 762-4000 or (+1) 203-925-4602

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.

