

## Click-iT® Plus EdU Imaging Kits

Search the handbook

Buy Now

Table 1 Contents and storage

Material	C10637	C10638	C10639	C10640	Concentration	Storage*
EdU (Component A)	5 mg	5 mg	5 mg	5 mg	NA	<ul style="list-style-type: none"> <li>• 2–8°C</li> <li>• Desiccate</li> <li>• Protect from light</li> <li>• DO NOT FREEZE</li> </ul>
Alexa Fluor® picolyl azide (Component B)	1 vial of Alexa Fluor® 488 picolyl azide (70 µL)	1 vial of Alexa Fluor® 555 picolyl azide (70 µL)	1 vial of Alexa Fluor® 594 picolyl azide (70 µL)	1 vial of Alexa Fluor® 647 picolyl azide (70 µL)	DMSO solution	
Dimethylsulfoxide (DMSO) (Component C)	4 mL	4 mL	4 mL	4 mL	NA	
Click-iT® EdU reaction buffer (Component D)	4 mL	4 mL	4 mL	4 mL	10X solution containing Tris-buffered saline	
Copper protectant (Component E)	1 vial, 0.5 mL	1 vial, 0.5 mL	1 vial, 0.5 mL	1 vial, 0.5 mL	NA	
Click-iT® EdU buffer additive (Component F)	400 mg	400 mg	400 mg	400 mg	NA	
Hoechst® 33342 (Component G)	35 µL	35 µL	35 µL	35 µL	10 mg/mL in water	

\*These storage conditions are appropriate when storing the entire kit upon receipt. For optimal storage conditions for each component, see vial labels. When stored as directed the product is stable for 1 year.

NA = Not applicable.

**Number of assays:** Sufficient material is supplied for 50 coverslips based on the protocol below.

**Approximate fluorescence excitation and emission maxima, in nm:** Alexa Fluor® 488 picolyl azide: 495/519; Alexa Fluor® 555 picolyl azide: 555/565; Alexa Fluor® 594 picolyl azide: 590/615; Alexa Fluor® 647 picolyl azide: 650/670; Hoechst® 33342: 350/461, bound to DNA.

## Introduction

Measuring a cell's ability to proliferate is a fundamental method for assessing cell health, determining genotoxicity, and evaluating anti-cancer drugs. The most accurate method is by directly measuring DNA synthesis. Initially, this was performed by incorporation of radioactive nucleosides (for example, <sup>3</sup>H-thymidine). This method was replaced by antibody-based detection of the nucleoside analog bromo-deoxyuridine (BrdU). The Click-iT® Plus EdU Assay is a novel alternative to the BrdU assay. EdU (5-ethynyl-2'-deoxyuridine) provided in the kit is a nucleoside analog of thymidine and is incorporated into DNA during active DNA synthesis<sup>1</sup>. Detection is based on a click reaction,<sup>2-5</sup> a copper catalyzed covalent reaction between a picolyl azide and an alkyne. In this application, the EdU contains the alkyne and the Alexa Fluor® dye contains the

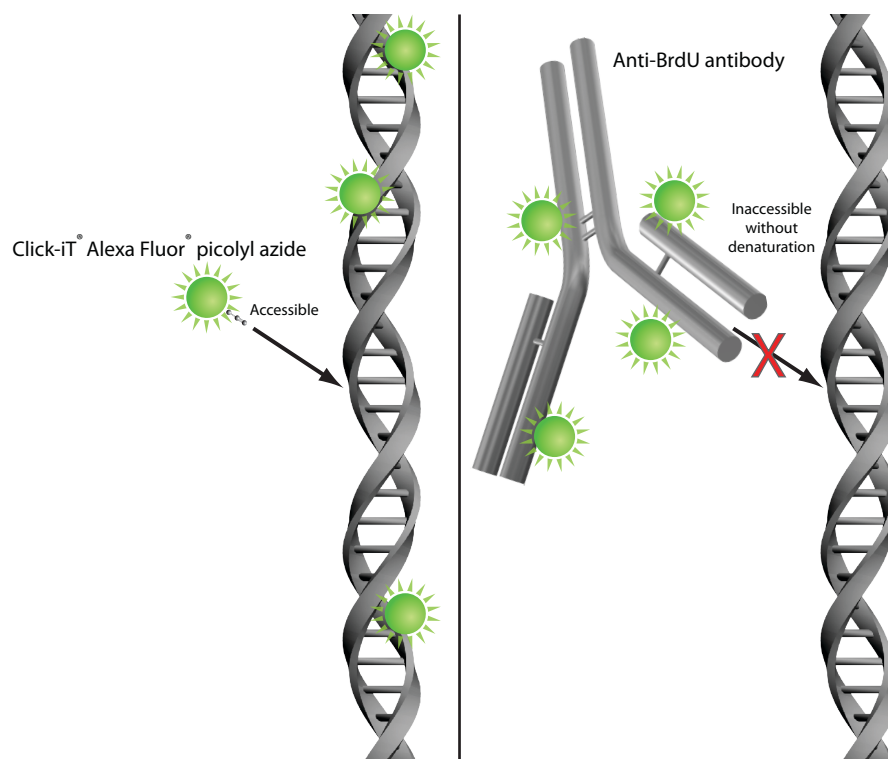
For Research Use Only. Not for use in diagnostic procedures.

picolyl azide. The advantages of the Click-iT<sup>®</sup> Plus EdU labeling are readily evident while performing the assay. The small size of the dye picolyl azide allows for efficient detection of the incorporated EdU using mild conditions. Standard aldehyde-based fixation and detergent permeabilization are sufficient for the Click-iT<sup>®</sup> Plus detection reagent to gain access to the DNA. This is in contrast to BrdU assays that require DNA denaturation (typically using HCl, heat, or digestion with DNase) to expose the BrdU so that it may be detected with an anti-BrdU antibody (Figure 1, below).

The denaturation step in the BrdU protocol can disrupt dsDNA integrity (which can affect nuclear counterstaining), and can also destroy cell morphology and antigen recognition sites. In contrast, the EdU assay kit is easy to use and fully compatible with DNA staining, including dyes for cell cycle analysis. This EdU assay kit can also be multiplexed with fluorescent proteins (for example, GFP, RFP, and mCherry; see Table 2 on page 3 for details). Unlike the BrdU assay, which relies upon antibodies which can exhibit nonspecific binding, the Click-iT<sup>®</sup> Plus EdU assay utilizes bioorthogonal (biologically unique) moieties, producing low backgrounds and high detection sensitivities.

The kit contains all of the components needed to label and detect the incorporated EdU, as well as perform cell cycle analysis on samples from adherent cells (Figure 2, page 3). For cell cycle analysis, the kit is supplied with blue fluorescent Hoechst<sup>®</sup> 33342 dye. The kits include sufficient reagents for labeling 50, 18 × 18 coverslips using 500 μL reaction buffer per test.

**Figure 1** Detection of the incorporated EdU with the Alexa Fluor<sup>®</sup> picolyl azide versus incorporated BrdU with an anti-BrdU antibody. The small size of the Alexa Fluor<sup>®</sup> picolyl azide eliminates the need to denature the DNA for the EdU detection reagent to gain access to the nucleotide.

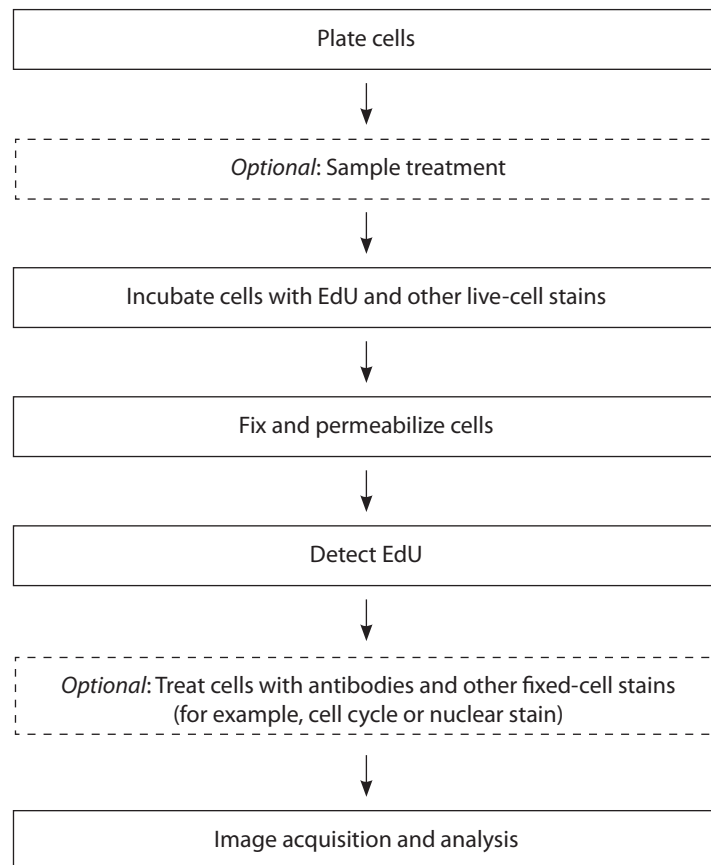


**Table 2** Click-iT® Plus detection reagent compatibility.

Molecule	Compatibility*
Qdot® nanocrystals	Use Qdot® nanocrystals <b>after</b> the Click-iT® Plus detection reaction.
Fluorescent proteins such as Green Fluorescent Protein (GFP)	<b>Completely compatible</b> with the Click-iT® Plus detection reaction.
Organic dyes such as Alexa Fluor® dyes, fluorescein (FITC)	<b>Completely compatible</b> with the Click-iT® Plus detection reaction.
Phalloidin	Phalloidin staining is <b>not compatible</b> with the Click-iT® Plus detection reaction. Use antibodies against other proteins, such as anti- $\alpha$ -tubulin, for visualization of the cytoskeleton.
Horseradish peroxidase (HRP)	Use HRP <b>after</b> the Click-iT® Plus detection reaction.
R-phycoerythrin (R-PE) and R-PE-tandems such as Alexa Fluor® 680-R-PE	<b>Completely compatible</b> with the Click-iT® Plus detection reaction.
Allophycocyanin (APC) and APC-tandems	<b>Completely compatible</b> with the Click-iT® Plus detection reaction.

\*Compatibility indicates whether the fluorescent molecule itself or the detection method involves components that are unstable in the presence of copper catalyst and copper protectant used for the Click-iT® Plus detection reaction.

**Figure 2** Workflow diagram for the Click-iT® Plus EdU Imaging Assay.



## Before You Begin

---

### Materials required but not provided

- Phosphate buffered saline (PBS, pH 7.2–7.6)
- Fixative (i.e., 3.7% Formaldehyde in PBS)
- Permeabilization reagent (i.e., 0.5% Triton<sup>®</sup> X-100 in PBS)
- 3% Bovine serum albumin (BSA) in PBS (3% BSA in PBS), pH 7.4
- Deionized water
- 18 × 18-mm coverslips
- Optional: 6-well microplate

### Cautions

- Hoechst<sup>®</sup> 33342 (Component G) is a known mutagen. Use the dye with appropriate precautions.
- DMSO (in Components B and C), provided as a solvent in this kit, is known to facilitate the entry of organic molecules into tissues. Handle reagents containing DMSO using equipment and practices appropriate for the hazards posed by such materials. Dispose of the reagents in compliance with all pertaining local regulations.

### Prepare stock solutions

1.1 Allow the vials to warm to room temperature before opening.

1.2 Prepare a 10 mM solution of EdU (Component A) by adding 2 mL DMSO (Component C) or an aqueous solution (i.e., buffer, saline) to EdU (Component A). Mix well.

After use, store any remaining stock solution at  $\leq -20^{\circ}\text{C}$ . When stored as directed the stock solution is stable for up to 1 year.

1.3 Prepare a working solution of 1X Click-iT<sup>®</sup> EdU reaction buffer (Component D): Transfer the solution (4 mL) in the Component D bottle to 36 mL of deionized water.

To make smaller amounts of 1X Click-iT<sup>®</sup> EdU reaction buffer, dilute volumes from the Component D bottle 1:10 with deionized water. After use, store any remaining 1X solution at 2–8°C. When stored as directed, this 1X solution is stable for 6 months.

1.4 To make a 10X stock solution of the Click-iT<sup>®</sup> EdU buffer additive (Component F): Add 2 mL deionized water to the vial, then mix until fully dissolved. After use, store any remaining stock solution at  $\leq -20^{\circ}\text{C}$ .

When stored as directed, this stock solution is stable for up to 1 year. If the solution develops a brown color, it has degraded and should be discarded.

### Label Cells with EdU

The following protocol was developed with A549, HeLa, and NIH/3T3 cells with an optimized EdU concentration of 10  $\mu$ M, but it can be adapted for any adherent cell type. Growth medium, cell density, cell type variations, and other factors may influence labeling. In initial experiments, we recommend testing a range of EdU concentrations to determine the optimal concentration for your cell type and experimental conditions. Although sufficient material is included with the kit for standard dose response, additional EdU (Cat. nos. A10044, E10187) is available. If currently using a BrdU based assay for cell proliferation, a similar concentration to BrdU is a good starting concentration for EdU.

- 2.1 Plate the cells on coverslips at the desired density and allow them to recover overnight before additional treatment.
- 2.2 Prepare a 2X working solution of EdU (Component A) in complete medium from the 10 mM stock solution. A suggested starting concentration is 10  $\mu$ M.
- 2.3 Prewarm the 2X EdU solution, then add an equal volume of the 2X EdU solution to the volume of media containing cells to be treated to obtain a 1X EdU solution (e.g., for a final concentration of 10  $\mu$ M, replace half of the media with fresh media containing 20  $\mu$ M EdU). We do not recommend replacing all of the media, because this could affect the rate of cell proliferation.
- 2.4 Incubate the cells for the desired length of time under conditions optimal for the cell type. The time of EdU exposure to the cells allows for direct measurement of cells synthesizing DNA. The choice of time points and the length of time depends on the cell growth rate. Pulse labeling of cells by brief exposures to EdU permits studies of cell-cycle kinetics.
- 2.5 Proceed immediately to **Fix and permeabilize cells** (steps 3.1–3.3, below) followed by **Detect EdU** (steps 4.1–4.7, page 6).

### Fix and permeabilize cells

**Note:** This protocol is optimized with a fixation step using 3.7% formaldehyde in PBS followed by a 0.5% Triton<sup>®</sup> X-100 permeabilization step, but it is also amenable to other fixation/permeabilization reagents such as methanol and saponin.

Transfer the coverslips into a 6-well plate for convenient processing such that each well contains a single coverslip.

- 3.1 After incubation, remove the media and add 1 mL of 3.7% formaldehyde in PBS to each well containing the coverslips. Incubate for 15 minutes at room temperature.
- 3.2 Remove the fixative and wash the cells in each well twice with 1 mL of 3% BSA in PBS.
- 3.3 Remove the wash solution. Add 1 mL of 0.5% Triton<sup>®</sup> X-100 in PBS to each well and incubate for 20 minutes at room temperature.

**Detect EdU** **Note:** This protocol uses 500  $\mu$ L of Click-iT<sup>®</sup> Plus reaction cocktail per coverslip. A smaller volume can be used as long as the remaining reaction components are maintained at the same ratios.

- 4.1 Prepare 1X Click-iT<sup>®</sup> EdU buffer additive (see Table 3, below) by diluting the 10X solution (prepared in step 1.4) 1:10 in deionized water. Prepare this solution **fresh** and use the solution on the same day.
- 4.2 Prepare Click-iT<sup>®</sup> Plus reaction cocktail according to Table 3 below. It is important to add the ingredients in the order listed in the table; otherwise, the reaction will not proceed optimally.

**Note:** Use the Click-iT<sup>®</sup> Plus reaction cocktail within 15 minutes of preparation.

**Table 3** Click-iT<sup>®</sup> Plus reaction cocktails.

Reaction components*	Number of coverslips						
	1	2	4	5	10	25	50
1X Click-iT <sup>®</sup> reaction buffer (prepared in step 1.3)	440 $\mu$ L	880 $\mu$ L	1.84 mL	2.25 mL	4.4 mL	10.9 mL	21.9 mL
Copper protectant (Component E)	10 $\mu$ L	20 $\mu$ L	40 $\mu$ L	50 $\mu$ L	100 $\mu$ L	250 $\mu$ L	500 $\mu$ L
Alexa Fluor <sup>®</sup> picolyl azide (Component B)	1.2 $\mu$ L	2.5 $\mu$ L	5 $\mu$ L	6 $\mu$ L	12.5 $\mu$ L	31 $\mu$ L	62 $\mu$ L
Reaction buffer additive (prepared in step 4.1)	50 $\mu$ L	100 $\mu$ L	200 $\mu$ L	250 $\mu$ L	500 $\mu$ L	1.25 mL	2.5 mL
Total volume	500 $\mu$ L	1 mL	2 mL	2.5 mL	5 mL	12.5 mL	25 mL

**\*Note:** Add the ingredients in the order listed in the table.

- 4.3 Remove the permeabilization buffer (step 3.3) and wash the cells in each well twice with 1 mL of 3% BSA in PBS. Remove the wash solution.
- 4.4 Add 0.5 mL of Click-iT<sup>®</sup> Plus reaction cocktail to each well containing a coverslip. Rock the plate briefly to insure that the reaction cocktail is distributed evenly over the coverslip.
- 4.5 Incubate the plate for 30 minutes at room temperature, **protected from light**.
- 4.6 Remove the reaction cocktail and wash each well once with 1 mL of 3% BSA in PBS. Remove the wash solution.

For nuclear staining, proceed to **Stain DNA** below. If no additional staining is desired, proceed to **Image and Analyze** on page 7.

- 4.7 *Optional:* Perform antibody labeling of the samples at this time, following the recommendations from the manufacturer of the primary and secondary antibody. It is important to keep the samples **protected from light** during incubations.

## Stain DNA

- 5.1 Wash each well with 1 mL of PBS. Remove the wash solution.
- 5.2 Dilute the Hoechst<sup>®</sup> 33342 (Component G) solution 1:2,000 in PBS to obtain a 1X Hoechst<sup>®</sup> 33342 solution (final concentration is 5  $\mu$ g/mL).

**Note:** A range between 2–10  $\mu$ g/mL of Hoechst<sup>®</sup> 33342 has been shown to work.

**5.3** Add 1 mL of 1X Hoechst® 33342 solution per well. Incubate for 30 minutes at room temperature, **protected from light**. Remove the Hoechst® 33342 solution.

**5.4** Wash each well twice with 1 mL of PBS. Remove the wash solution.

**Image and analyze** Click-iT® Plus EdU cells are compatible with all methods of slide preparation including wet mount or prepared mounting media. See Table 4 below for the approximate fluorescence excitation/emission maxima for Alexa Fluor® dyes and Hoechst® 33342 dye bound to DNA.

**Table 4** Approximate fluorescence excitation/emission maxima.

Fluorophore	Excitation (nm)	Emission (nm)
Alexa Fluor® 488 picolyl azide	495	519
Alexa Fluor® 555 picolyl azide	555	565
Alexa Fluor® 594 picolyl azide	590	615
Alexa Fluor® 647 picolyl azide	650	670
Hoechst® 33342, bound to DNA	350	461

## References

**1.** Proc Natl Acad Sci USA 105, 2415 (2008); **2.** ChemBioChem 4, 1147 (2003); **3.** J Am Chem Soc 125, 3192 (2003); **4.** Angew Chem Int Ed Engl 41, 2596 (2002); **5.** Angew Chem Int Ed Engl 40, 2004 (2001).

## Product List

Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
<a href="#">Buy Now</a> C10637	Click-iT® Plus EdU Alexa Fluor® 488 Imaging Kit *for 50 coverslips*	1 kit
<a href="#">Buy Now</a> C10638	Click-iT® Plus EdU Alexa Fluor® 555 Imaging Kit *for 50 coverslips*	1 kit
<a href="#">Buy Now</a> C10639	Click-iT® Plus EdU Alexa Fluor® 594 Imaging Kit *for 50 coverslips*	1 kit
<a href="#">Buy Now</a> C10640	Click-iT® Plus EdU Alexa Fluor® 647 Imaging Kit *for 50 coverslips*	1 kit
<b>Related Products</b>		
<a href="#">Buy Now</a> C10289	Click-iT® AHA Alexa Fluor® 488 Protein Synthesis HCS Assay	1 kit
<a href="#">Buy Now</a> C10429	Click-iT® HPG Alexa Fluor® 594 Protein Synthesis Assay Kit	1 kit
<a href="#">Buy Now</a> C10428	Click-iT® HPG Alexa Fluor® 488 Protein Synthesis Assay Kit	1 kit
<a href="#">Buy Now</a> C10327	Click-iT® RNA Alexa Fluor® 488 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10328	Click-iT® RNA Alexa Fluor® 594 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10329	Click-iT® RNA Alexa Fluor® 488 Imaging Kit *for 25 coverslips*	1 kit
<a href="#">Buy Now</a> C10330	Click-iT® RNA Alexa Fluor® 594 Imaging Kit *for 25 coverslips*	1 kit
<a href="#">Buy Now</a> C10350	Click-iT® EdU Alexa Fluor® 488 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10351	Click-iT® EdU Alexa Fluor® 488 HCS Assay *10-plate size*	1 kit
<a href="#">Buy Now</a> C10352	Click-iT® EdU Alexa Fluor® 555 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10353	Click-iT® EdU Alexa Fluor® 555 HCS Assay *10-plate size*	1 kit
<a href="#">Buy Now</a> C10354	Click-iT® EdU Alexa Fluor® 594 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10355	Click-iT® EdU Alexa Fluor® 594 HCS Assay *10-plate size*	1 kit
<a href="#">Buy Now</a> C10356	Click-iT® EdU Alexa Fluor® 647 HCS Assay *2-plate size*	1 kit
<a href="#">Buy Now</a> C10357	Click-iT® EdU Alexa Fluor® 647 HCS Assay *10-plate size*	1 kit
<a href="#">Buy Now</a> A10044	EdU [5-ethynyl-2'-deoxyuridine]	50 mg
<a href="#">Buy Now</a> E10187	EdU [5-ethynyl-2'-deoxyuridine]	500 mg
<a href="#">Buy Now</a> E10415	EdU [5-ethynyl-2'-deoxyuridine]	5 g
<a href="#">Buy Now</a> H3570	Hoechst® 33342, trihydrochloride, trihydrate *10 mg/mL solution in water*	10 mL

# Purchaser Notification

---

## Corporate Headquarters

5791 Van Allen Way  
Carlsbad, CA 92008  
USA  
Phone: +1 760 603 7200  
Fax: +1 760 602 6500  
Email: [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

## European Headquarters

Inchinnan Business Park  
3 Fountain Drive  
Paisley PA4 9RF  
UK  
Phone: +44 141 814 6100  
Toll-Free Phone: 0800 269 210  
Toll-Free Tech: 0800 838 380  
Fax: +44 141 814 6260  
Tech Fax: +44 141 814 6117  
Email: [euroinfo@invitrogen.com](mailto:euroinfo@invitrogen.com)  
Email Tech: [eurotech@invitrogen.com](mailto:eurotech@invitrogen.com)

## Japanese Headquarters

LOOP-X Bldg. 6F  
3-9-15, Kaigan  
Minato-ku, Tokyo 108-0022  
Japan  
Phone: +81 3 5730 6509  
Fax: +81 3 5730 6519  
Email: [jpinfo@invitrogen.com](mailto:jpinfo@invitrogen.com)

Additional international offices are listed at  
[www.lifetechnologies.com](http://www.lifetechnologies.com)

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

## Obtaining Support

For the latest services and support information for all locations, go to [www.lifetechnologies.com](http://www.lifetechnologies.com).

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support ([techsupport@lifetech.com](mailto:techsupport@lifetech.com))
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

## SDS

Safety Data Sheets (SDSs) are available at [www.lifetechnologies.com/sds](http://www.lifetechnologies.com/sds).

## Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

## Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

## Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

## Important Licensing Information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners. Triton is a registered trademark of Union Carbide Corporation. Hoechst is a registered trademark of Hoechst GmbH.

©2014 Life Technologies Corporation. All rights reserved.