


Click-iT™ Plus EdU Flow Cytometry Assay Kits

Catalog Numbers C10632, C10634, C10636, C10645, and C10646

Pub. No. MAN0009883 **Rev.** E.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

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, i.e., ³H-thymidine. This method was replaced by antibody-based detection of the nucleoside analog bromo-deoxyuridine (BrdU). The Click-iT™ Plus EdU Flow Cytometry Assay Kits are novel alternatives to the BrdU assay. EdU (5-ethynyl-2'-deoxyuridine) is a nucleoside analog to thymidine and is incorporated into DNA during active DNA synthesis. Detection is based on a click reaction, a copper catalyzed covalent reaction between a picolyl azide and an alkyne. In this application, the alkyne is found in the ethynyl moiety of EdU, while the picolyl azide is coupled to Alexa Fluor™ 350 dye, Alexa Fluor™ 488 dye, Alexa Fluor™ 594 dye, Alexa Fluor™ 647 dye, or Pacific Blue™ dye. Standard flow cytometry methods are used for determining the percentage of S-phase cells in the population (see "Expected results" on page 5).

The advantage of Click-iT™ Plus EdU labeling is that the small size of the picolyl dye 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™ Plus detection reagent to gain access to the DNA. This is in contrast to BrdU assays that require DNA denaturation (using acid, heat, or digestion with DNase) to expose the BrdU so that it may be detected with an anti-BrdU antibody. Sample processing for the BrdU assay can result in signal alteration of the cell cycle distribution as well as the destruction of antigen recognition sites when using the acid denaturation method. In contrast, the Click-iT™ Plus EdU cell proliferation kit is compatible with cell cycle dyes, R-PE, R-PE tandems, and fluorescent proteins such as GFP, RFP, and mCherry. See page 5 for examples of Click-iT™ Plus EdU Alexa Fluor™ 488 multiplexed with R-PE, Click-iT™ Plus EdU Alexa Fluor™ 350 multiplexed with Pacific Orange™ and FxCycle™ Far Red, and Click-iT™ Plus EdU Alexa Fluor™ 647 multiplexed with GFP and FxCycle™ Violet. The EdU assay can also be multiplexed with antibodies against surface and intracellular markers (see Table 3).

Contents and storage

Upon receipt, store the kit at 2°C to 8°C. For optimal storage conditions for each component, see vial labels. When stored as directed the product is stable for up to 1 year after receipt.

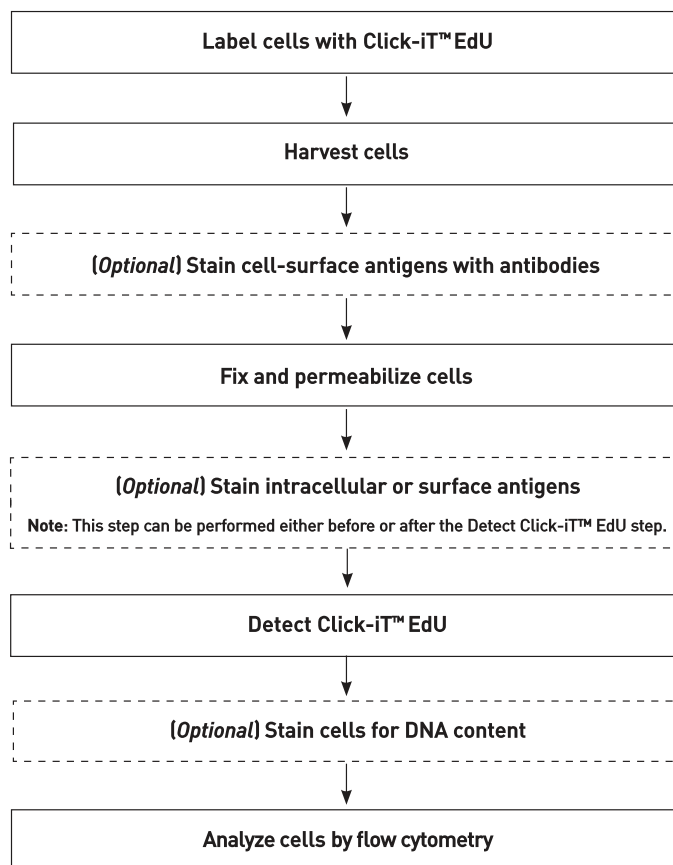
Contents	C10632	C10634	C10636	C10645	C10646	Storage
EdU (Component A)	10 mg					<ul style="list-style-type: none"> • 2–8°C • Desiccate • Protect from light • Do not freeze
Alexa Fluor™ 350 picolyl azide; in DMSO (Component B)	—	—	—	130 µL	—	
Alexa Fluor™ 488 picolyl azide; in DMSO (Component B)	130 µL	—	—	—	—	
Alexa Fluor™ 594 picolyl azide; in DMSO (Component B)	—	—	—	—	130 µL	
Alexa Fluor™ 647 picolyl azide; in DMSO (Component B)	—	130 µL	—	—	—	
Pacific Blue™ picolyl azide; in DMSO (Component B)	—	—	130 µL	—	—	
Dimethylsulfoxide [DMSO] (Component C)	4.5 mL					
Click-iT™ fixative [4% paraformaldehyde in PBS] (Component D)	5 mL					
Click-iT™ fixative saponin-based permeabilization and wash reagent, 10X (Component E)	50 mL					
Copper protectant, 100 mM aqueous solution (Component F)	0.5 mL					
Click-iT™ EdU buffer additive (Component G)	400 mg					

Required materials not supplied

- 1% bovine serum albumin (BSA) in phosphate buffered saline (PBS), pH 7.1–pH 7.4
- Buffered saline solution, such as PBS, D-PBS, or TBS
- Deionized water or 18 MΩ purified water
- 12 × 75 mm round bottom tubes, or other flow cytometry tubes

Workflow

The following protocol was developed with Jurkat cells, a human T cell line, and using an EdU concentration of 10 μ M, and can be adapted for any 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. If currently using a BrdU based assay for cell proliferation, a similar concentration to BrdU is a good starting concentration for EdU. If using whole blood as the sample, we recommend heparin as the anticoagulant for collection.



Procedural guidelines



CAUTION!

- 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.
- Click-iT™ fixative (Component D) contains paraformaldehyde, which is harmful. Use with appropriate precautions.
- Click-iT™ saponin-based permeabilization and wash reagent (Component E) contains sodium azide, which yields highly toxic hydrazoic acid under acidic conditions. Dilute azide compounds in running water before discarding to avoid accumulation of potentially explosive deposits in plumbing.

Allow vials to warm to room temperature before opening them to prepare reagents.

Prepare 10 mM EdU solution

Add 4 mL of DMSO (Component C) or aqueous solution (PBS) to Component A and mix well to prepare a 10 mM solution of EdU. 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.

Prepare 10X Click-iT™ EdU buffer additive

Add 2 mL of deionized water to the vial and mix until the Click-iT™ EdU buffer additive is fully dissolved. 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.

Prepare 1X Click-iT™ permeabilization and wash reagent

Note: There is a degree of variability that has been observed in the color of the saponin solution (component E), from light yellow to dark brown. There is no difference in the kit performance by flow cytometry for the different colors of saponin that are observed.

add 50 mL of Component E to 450 mL of 1% BSA in PBS to prepare 500 mL of 1X Click-iT™ saponin-based permeabilization and wash reagent.

Smaller amounts can be prepared by diluting a volume of Component E 1:10 with 1% BSA in PBS. After use, store any remaining solutions at 2–8°C. When stored as directed, the 1X solution is stable for 6 months and the 10X solution is stable for 12 months after receipt.

Label cells with Click-iT™ Edu

1. Suspend the cells in an appropriate tissue culture medium to obtain optimal conditions for cell growth. Disturbing the cells by temperature changes or washing prior to incubation with EdU slows the growth of the cells during incorporation.
2. Add EdU to the culture medium at the desired final concentration and mix well. We recommend a starting concentration of 10 μM for 1–2 hours. For longer incubations, use lower concentrations. For shorter incubations, higher concentrations may be required. For a negative staining control, include cells from the same population that have not been treated with EdU.
3. Incubate under conditions optimal for cell type for the desired length of time. Altering the amount of time the cells are exposed to EdU or subjecting the cells to pulse labeling with EdU allows the evaluation of various DNA synthesis and proliferation parameters. Effective time intervals for pulse labeling and the length of each pulse depend on the cell growth rate.
4. Harvest cells and proceed immediately to "Stain cell-surface antigens with antibodies" if performing antibody surface labeling; otherwise continue to "Fix and permeabilize cells".

(Optional) Stain cell-surface antigens with antibodies

1. Wash cells once with 3 mL of 1% BSA in PBS, pellet cells by centrifugation, and remove supernatant.
2. Dislodge the pellet and resuspend cells at 1×10^7 cells/mL in 1% BSA in PBS.
3. Add 100 μL of cell suspension or whole blood sample to flow tubes.
4. Add surface antibodies and mix well (see "Performance characteristics" on page 5).
Note: Do not use Qdot™ antibody conjugates before performing the click reaction; see "Stain intracellular or surface antigens" to label with Qdot™ fluorophores.
5. Incubate for the recommended time and temperature, protected from light.
6. Proceed to "Fix and permeabilize cells".

Fix and permeabilize cells

The Click-iT™ saponin-based permeabilization and wash reagent can be used with whole blood or cell suspensions containing red blood cells, as well as with cell suspensions containing more than one cell type. This permeabilization and wash reagent maintains the morphological light scatter characteristics of leukocytes while lysing red blood cells.

1. Wash the cells once with 3 mL of 1% BSA in PBS, pellet the cells, and remove the supernatant.
2. Dislodge the pellet, add 100 μL of Click-iT™ fixative (Component D), and mix well.
3. Incubate the cells for 15 minutes at room temperature, protected from light.
4. Wash the cells with 3 mL of 1% BSA in PBS, pellet the cells, and remove the supernatant. Repeat the wash step if red blood cells or hemoglobin are present in the sample. Remove all residual red blood cell debris and hemoglobin before proceeding.
5. Dislodge the cell pellet and resuspend the cells in 100 μL of 1X Click-iT™ permeabilization and wash reagent (see page 2), and mix well. Incubate the cells for 15 minutes then proceed directly to "Detect Click-iT™ Edu" on page 4, or proceed to "Stain intracellular or surface antigens".

(Optional) Stain intracellular or surface antigens

This step can be performed either before or after the Detect Click-iT™ Edu step.

1. Add antibodies against intracellular antigens or against surface antigens that use Qdot™ antibody conjugates. Mix well.
2. Incubate the tubes for the time and temperature required for antibody staining, protected from light.
3. Wash the cells according to the following conditions, depending on whether you have already performed the Detect Click-iT™ Edu step.

Click-iT™ Edu detection	Action
Already performed the Detect Click-iT™ Edu step.	<ol style="list-style-type: none">1. Wash each tube with 3 mL of 1X Click-iT™ permeabilization and wash reagent.2. Centrifuge the cells, and remove the supernatant.3. Dislodge the cell pellet and resuspend the cells in 500 μL of 1X Click-iT™ permeabilization and wash reagent.4. Proceed to "Stain cells for DNA content" or "Analyze cells by flow cytometry" on page 4.
Have not yet performed the Detect Click-iT™ Edu step.	<ol style="list-style-type: none">1. Wash the cells once with 3 mL of 1% BSA in PBS, pellet the cells, and remove the supernatant.2. Dislodge the cell pellet and add 100 μL of 1X Click-iT™ permeabilization and wash reagent and mix well.3. Proceed to "Detect Click-iT™ Edu" on page 4.

Detect Click-iT™ EdU

1. Prepare 1X Click-iT™ EdU buffer additive by diluting the 10X stock solution (see page 2) 1:10 in deionized water.
2. Prepare the Click-iT™ Plus reaction cocktail according to the following table.

Table 1

Reaction components	Number of reactions						
	1	2	5	10	15	30	50
PBS, D-PBS, or TBS	438 µL	875 µL	2.19 mL	4.38 mL	6.57 mL	13.2 mL	21.9 mL
Copper protectant (Component F)	10 µL	20 µL	50 µL	100 µL	150 µL	300 µL	500 µL
Fluorescent dye picolyl azide	2.5 µL	5 µL	12.5 µL	25 µL	37.5 µL	75 µL	125 µL
Reaction Buffer Additive (prepared in step 1)	50 µL	100 µL	250 µL	500 µL	750 µL	1.5 mL	2.5 mL
Total reaction volume	500 µL	1 mL	2.5 mL	5 mL	7.5 mL	15 mL	25 mL

Note: Use the Click-iT™ Plus reaction cocktail within 15 minutes of preparation.

3. Add 0.5 mL of Click-iT™ Plus reaction cocktail to each tube of fixed cells and mix well.

Note: The total volume for each reaction mixture is 600 µL.

4. Incubate the reaction mixture for 30 minutes at room temperature, protected from light.
5. Wash the cells once with 3 mL of 1X Click-iT™ permeabilization and wash reagent (see page 2), centrifuge the cells, and remove the supernatant.

Option	Description
If proceeding to "Analyze cells by flow cytometry"	Go to page 4.
If proceeding to "Stain intracellular or surface antigens"	Dislodge the cell pellet and resuspend the cells in 100 µL of 1X Click-iT™ permeabilization and wash reagent.
If proceeding to "Stain cells for DNA content"	Add 500 µL of 1X Click-iT™ permeabilization and wash reagent.

(Optional) Stain cells for DNA content

1. If necessary, add Ribonuclease A to each tube and mix (see the following table).
2. Add the appropriate DNA stain to each tube, mix well, then incubate as recommended for each DNA stain.

Table 2

DNA content stain	RNase required?	Click-iT™ Plus EdU stain compatibility				
		Alexa Fluor™ 350 picolyl azide	Alexa Fluor™ 488 picolyl azide	Alexa Fluor™ 594 picolyl azide	Alexa Fluor™ 647 picolyl azide	Pacific Blue™ picolyl azide
FxCycle™ PI/RNase	No	Yes	Yes	No	Yes	Yes
FxCycle™ Violet	No	No	Yes	Yes	Yes	No
FxCycle™ Far Red	Yes	Yes	Yes	Yes	No	Yes
SYTOX™ AADvanced™ stain	Yes	Yes	Yes	Yes	Yes	Yes
Propidium iodide (PI)	Yes	Yes	Yes	No	Yes	Yes

Analyze cells by flow cytometry

If measuring total DNA content on a traditional flow cytometer using hydrodynamic focusing, use a low flow rate during acquisition. If using the Attune™ or Attune™ NxT Flow Cytometer, all collection rates may be used without loss of signal integrity if the event rate is kept below 10,000 events per second. However, for each sample within an experiment, the same collection rate and cell concentration should be used. The fluorescent signal generated by DNA content stains is best detected with linear amplification. The fluorescent signal generated by Click-iT™ Plus EdU labeling is best detected with logarithmic amplification.

Analyze the cells using a flow cytometer.

For detection of EdU with	Use
Alexa Fluor™ 350 picolyl azide	350 nm excitation with a 440 emission filter (450/50 nm or similar)
Alexa Fluor™ 488 picolyl azide	488 nm excitation with a green emission filter (530/30 nm or similar)
Alexa Fluor™ 594 picolyl azide	532 or 561 nm excitation with a 620 emission filter (620/15 nm or similar)
Alexa Fluor™ 647 picolyl azide	633/635 nm excitation with a red emission filter (660/20 nm or similar)
Pacific Blue™ picolyl azide	405 nm excitation with a violet emission filter (450/40 nm or similar)

Performance characteristics

The EdU assay can also be multiplexed with antibodies against surface and intracellular markers

Compatibility indicates whether the fluorescent molecule itself or the detection methods involve components that are unstable in the presence of the copper catalyst used for the Click-iT™ Plus EdU detection reaction.

Table 3 Click-iT™ Plus EdU detection reagent compatibility

Fluorescent molecule	Compatibility
R-phycoerythrin (R-PE) and R-PE based tandems (i.e., Alexa Fluor™ 610-RPE)	Compatible, except Alexa Fluor™ 594 with R-PE-Cy7™ tandems
Fluorescent proteins (GFP)	Compatible
PerCP, allophycocyanin (APC) and APC-based tandems (i.e., Alexa Fluor™ 680-APC)	Compatible
Organic dyes such as Alexa Fluor™ dyes, fluorescein (FITC)	Compatible
Qdot™ nanocrystals	Use Qdot™ nanocrystals after the Click-iT™ Plus detection reaction.

Expected results

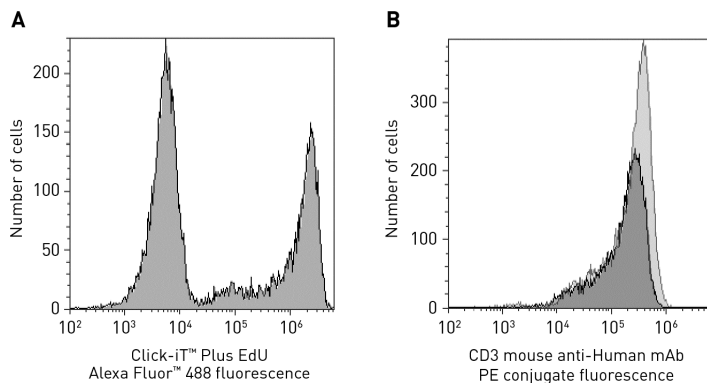


Fig. 1 Fluorescence signal from Alexa Fluor™ 488 Click-iT™ Plus EdU Flow Cytometry Assay Kits and CD3 mouse anti-Human mAb PE conjugate

Jurkat (human T-cell leukemia) cells were treated with 10 μ M EdU for 2 hours, stained with CD3 mouse anti-Human mAb PE conjugate (Cat. no. MHCD0304) and detected according to the recommended staining protocol. The figures show a clear separation of proliferating cells which have incorporated EdU and nonproliferating cells which have not. Panel A shows data from cells labeled with Alexa Fluor™ 488 picolyl azide analyzed on an Attune™ Acoustic Focusing Cytometer using 488 nm excitation and a 530/30 nm bandpass emission filter; Panel B shows the same cells using 635 nm excitation and a 574/26 nm bandpass emission filter. The black outlined histogram is the cells stained with CD3 mouse anti-Human mAb PE conjugate and Click-iT™ Plus EdU Alexa Fluor™ 488 picolyl azide. The gray outlined histogram is the CD3 mouse anti-Human mAb PE conjugate positive control cells treated the same but without copper in the reaction.

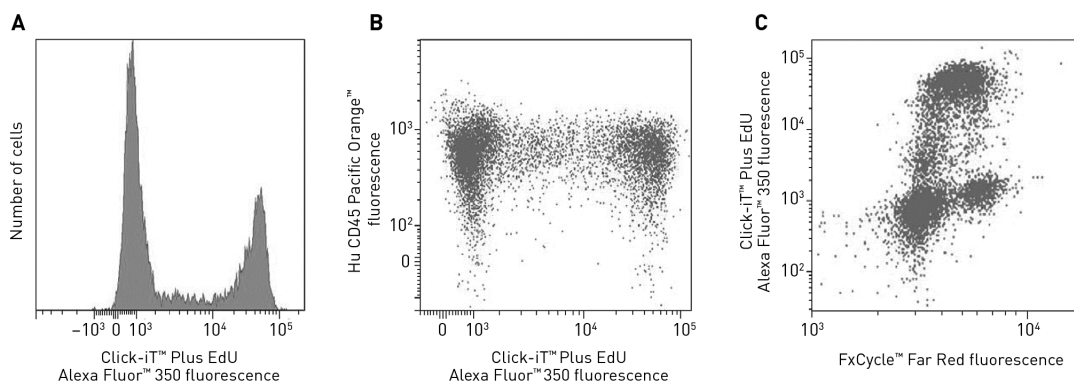


Fig. 2 Fluorescence signal and dual parameter plots from Alexa Fluor™ 350 Click-iT™ Plus EdU Flow Cytometry Assay Kit, CD45-Pacific Orange™ and FxCycle™ Far Red

Jurkat (human T-cell leukemia) cells were treated with 10 μ M EdU for 2 hours, stained with CD45-Pacific Orange™ (Cat. no. MHCD4530), and detected according to the recommended staining protocol. The figures show a clear separation of proliferating cells which have incorporated EdU and non-proliferating cells which have not. Panel A shows data from cells labeled with Alexa Fluor™ 350 picolyl azide analyzed on a BD™ LSRII flow cytometer using UV excitation and a 450/50 nm bandpass emission filter; Panel B shows the same cells using 488 nm excitation and a 530/30 nm bandpass emission filter for detection of the CD45-Pacific Orange™ and UV excitation, and a 450/50 nm bandpass emission filter for detection of the Alexa Fluor™ 350 picolyl azide; Panel C shows the dual parameter plot of the Click-iT™ Plus EdU Alexa Fluor™ 350 and FxCycle™ Far Red. Data were collected and analyzed using a BD™ LSRII flow cytometer using UV excitation and a 450/50 nm bandpass

emission filter for detection of the Alexa Fluor™ 350 picolyl azide and 633 nm excitation and a 660/20 nm bandpass emission filter for detection of the FxCycle™ Far Red fluorescence. This figure combines DNA content with EdU; cells that are positive for both labels are in S-phase of the cell cycle.

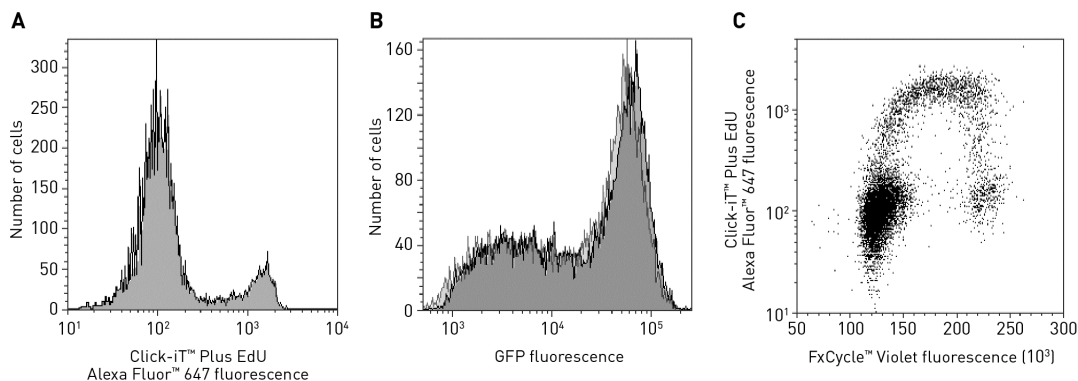


Fig. 3 Fluorescence signal and dual parameter plot of Alexa Fluor™ 647 Click-iT™ Plus EdU Flow Cytometry Assay Kits, Green Fluorescent Protein, and FxCycle™ Violet

A375 (human malignant melanoma) cells were treated with 10 µM EdU for 2 hours and detected according to the recommended staining protocol. The figures show a clear separation of proliferating cells which have incorporated EdU and nonproliferating cells which have not. Panel A shows data from cells labeled with Alexa Fluor™ 647 picolyl azide analyzed on an Attune™ Acoustic Focusing Cytometer using 635 nm excitation and a 660/20 nm bandpass emission filter; Panel B shows the same cells using 488 nm excitation and a 530/30 nm bandpass emission filter. The black outlined histogram is the GFP-expressing cells and Click-iT™ Plus EdU Alexa Fluor™ 647 picolyl azide. The gray outlined histogram is the GFP-expressing positive control cells treated the same but without copper in the reaction. Panel C shows the dual parameter plot of the Click-iT™ Plus EdU Alexa Fluor™ 647 and FxCycle™ Violet. Data were collected and analyzed using an Attune™ Acoustic Cytometer using 635 nm excitation and a 660/20 nm bandpass emission filter for detection of the EdU Alexa Fluor™ 647 picolyl azide and 405 nm excitation and a 450/40 bandpass emission filter for detection of the FxCycle™ Violet fluorescence. This figure combines DNA content with EdU; cells that are positive for both labels are in S-phase of the cell cycle.

Product list

Product	Amount	Cat. No.
Click-iT™ Plus EdU Alexa Fluor™ 350 Flow Cytometry Assay Kit	1 kit (50 assays)	C10645
Click-iT™ Plus EdU Alexa Fluor™ 488 Flow Cytometry Assay Kit	1 kit (50 assays)	C10632
Click-iT™ Plus EdU Alexa Fluor™ 594 Flow Cytometry Assay Kit	1 kit (50 assays)	C10646
Click-iT™ Plus EdU Alexa Fluor™ 647 Flow Cytometry Assay Kit	1 kit (50 assays)	C10634
Click-iT™ Plus EdU Pacific Blue™ Flow Cytometry Assay Kit	1 kit (50 assays)	C10636
Click-iT™ Plus EdU Alexa Fluor™ 647 Flow Cytometry Assay Kit	1 kit (100 assays)	C10635
Click-iT™ Plus EdU Alexa Fluor™ 488 Flow Cytometry Assay Kit	1 kit (100 assays)	C10633
EdU (5-ethynyl-2'-deoxyuridine)	50 mg	A10044
BrdU mouse monoclonal antibody (Clone MoBU-1) - Pacific Blue™, for flow cytometry	1 each (100 tests)	B35129
BrdU mouse monoclonal antibody (Clone MoBU-1) - Alexa Fluor™ 488, for flow cytometry	1 each (100 tests)	B35139
BrdU mouse monoclonal antibody (Clone MoBU-1) - Alexa Fluor™ 647, for flow cytometry	1 each (100 tests)	B35140
BrdU mouse monoclonal antibody (Clone MoBU-1) - unconjugated, for flow cytometry	1 each (100 tests)	B35141
Click-iT™ EdU Alexa Fluor™ 488 HCS Assay	1 kit (2 plates)	C10350
Click-iT™ EdU Alexa Fluor™ 488 HCS Assay	1 kit (10 plates)	C10351
Click-iT™ EdU Alexa Fluor™ 594 HCS Assay	1 kit (2 plates)	C10354
Click-iT™ EdU Alexa Fluor™ 647 HCS Assay	1 kit (2 plates)	C10356
FxCycle™ Violet Stain, for flow cytometry	1 kit (500 assays)	F10347
FxCycle™ Far Red Stain, for flow cytometry	1 kit (500 assays)	F10348
FxCycle™ PI/RNase Staining Solution	100 mL (200 assays)	F10797
Hoechst™ 33342, trihydrochloride, trihydrate (10 mg/mL solution in water)	10 mL	H3570
Propidium iodide - 1.0 mg/mL solution in water	10 mL	P3566
SYTOX™ AADvanced™ dead cell stain, for flow cytometry at 488 excitation	1 kit (100 tests)	S10349
Vybrant™ DyeCycle™ Violet stain (5 mM in water)	200 µL (200 assays)	V35003
RNase A (20 mg/mL)	10 mL	12091-021
Dulbecco's Phosphate Buffered Saline 1X without Calcium Chloride without Magnesium Chloride	500 mL	14190-144
	10 × 500 mL	14190-250

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.