
BrdU Staining Kit for Flow Cytometry

Research Use Only

Introduction

The eBioscience BrdU Staining Kit for Flow Cytometry contains the necessary reagents and buffers for identifying and examining proliferating cells by flow cytometric analysis. Cycling cells are incubated with 5-bromo-2'-deoxyuridine (BrdU), a synthetic analog of thymidine which incorporates into newly synthesized genomic DNA during the S-phase of mitosis. Following DNA denaturation, the cells are stained for BrdU incorporation along with any other cell surface and/or intracellular targets of interest.

General Notes

Please use caution when handling. Product contains formaldehyde.

Protocol

Materials Provided

- BrdU (32.5 mM, 10 mg/mL): 5 x 1 mL sterile vials; store at less than or equal to -70°C. Avoid multiple freeze-thaws. Vial should be opened under sterile conditions.
- DNase I (1 mg/mL; 0.3 mg/vial): 10 x 0.3 mL vials; store at less than or equal to -70°C. Each vial can be used to treat 10 samples. Avoid multiple freeze-thaws.
- Anti-BrdU Antibody (clone BU20A), fluorochrome-conjugated: 1 x 100 test vial; store at 2-8°C.
- BrdU Staining Buffer Concentrate (4X): 1 x 30 mL bottle; store at 2-8°C. This buffer contains formaldehyde. Please handle appropriately.
- Fixation/Permeabilization Diluent: 1 x 100 mL bottle; store at 2-8°C.

Other Materials Needed

- Sterile 1X PBS
- Flow Cytometry Staining Buffer (eBioscience Cat. No. 00-4222)
- 12 x 75 mm round bottom test tubes
- Optional:
 - Primary antibodies (directly conjugated)
 - Fixable Viability Dye (FVD) eFluor® 450 (eBioscience Cat. No. 65-0863), eFluor® 660 (eBioscience Cat. No. 65-0864), eFluor® 780 (eBioscience Cat. No. 65-0865), eFluor® 506 (eBioscience Cat. No. 65-0866), eFluor® 520 (eBioscience Cat. No. 65-0867) and eFluor® 455UV (eBioscience Cat. No. 65-0868)

Note: The antibodies used for surface staining can be added after BrdU and Fixable Viability Dye labeling (but before fixation). Alternatively if the antibody(s) is known to recognize a formaldehyde-fixed epitope, it can be added concurrently with the BrdU antibody.

BrdU Staining Buffer Working Solution Preparation

- Prepare fresh 1X BrdU Staining Buffer working solution by diluting BrdU Staining Buffer Concentrate (1 part) with Fixation/Permeabilization Diluent (3 parts). Mix by gentle inversion, do not vortex. You will need 1 mL of the 1X BrdU Staining Buffer working solution for each sample. Use caution and handle appropriately as the buffer contains formaldehyde.

For additional questions, please contact Technical Support at +1-888-810-6168 (US) or +43 1 796 4040 120 (Europe/International), or send us an email at Tech_Support@affymetrix.com

BrdU Staining Kit for Flow Cytometry

Research Use Only

Experimental Procedure

Step 1: In vitro labeling of 10^5 to 10^8 dividing cells with 10 μ M BrdU for 45 minutes at 37°C.

- Under sterile conditions thaw BrdU on ice and dilute to a working concentration of 1 mM with sterile 1X PBS.
- Add 10 μ M BrdU to each sample. (For example, add 10 μ L of 1 mM BrdU directly to every milliliter of tissue culture medium.)
- Incubate your cells long enough to allow incorporation of BrdU. The timing will be dependent on your culture conditions (e.g., stimulants used) and the proliferation kinetics of your cells. Therefore the incubation time will need to be determined empirically. After the incubation, harvest the cells.
- Wash cells by adding 2 mL of Flow Cytometry Staining Buffer (or azide-free PBS if proceeding to Step 2) and then centrifuge at 300-400 xg for 5 minutes at room temperature. Discard the supernatant.

Step 2: [Optional] Stain with Fixable Viability Dye (FVD) to label dead cells before fixation.

Note: Allow the vial of Fixable Viability Dye to equilibrate to room temperature before opening. The dye must be used with azide-free PBS. For consistent staining of cells in tubes, do not stain in less than 0.5 mL. Please refer to the eBioscience website Best Protocols “[Viability Staining Protocol, Protocol C: Staining Dead Cells with eBioscience Fixable Viability eFluor® Dyes](#)” for additional information. (Proceed to Step 3 if a FVD will not be used.)

- Wash cells one additional time with 2 mL of azide-free PBS, as described in Step 1d.
- Resuspend cells at $1-10 \times 10^6$ cells/mL in azide-free PBS.
- Add 1 μ L of Fixable Viability Dye per 1 mL of cells and vortex immediately.
- Incubate for 30 minutes at 2-8°C in the dark.
- Wash cells 1-2 times with Flow Cytometry Staining Buffer, as described in Step 1d.
- Resuspend cells at $1-10 \times 10^6$ cells/mL in Flow Cytometry Staining Buffer.

Step 3: [Optional] Stain cell surface antigen(s).

Note: For additional information, please refer to the eBioscience website Best Protocols “[Staining Cell Surface Antigens for Flow Cytometry](#).” (Proceed to Step 4 if cell surface antigens will not be examined or if the antibody(s) is known to recognize a formaldehyde-fixed epitope.)

- Aliquot 50 μ L of cell suspension to each tube or well. The cell number should be determined empirically but can range from 10^5 to 10^8 cells/test per tube.
- Add the recommended amount (refer to the Technical Data Sheet for each product) of each fluorochrome-conjugated primary antibody(s) in an appropriate volume of Flow Cytometry Staining Buffer such that the final staining volume is 100 μ L. (For example, add 50 μ L of an antibody mix to 50 μ L of cells.) Mix gently.
- Incubate for at least 30 minutes at 2-8°C in the dark.
- Wash the cells twice with Flow Cytometry Staining Buffer, as described in Step 1d.

Step 4: Fix cells and intracellularly stain with Anti-BrdU.

- Thaw DNase I solution on ice. Once thawed, prepare a working solution of DNase I by adding 300 μ L of the DNase I solution to 700 μ L of Flow Cytometry Staining Buffer and mix gently. Store on ice until ready for use in Step 4f, below.
- If cells were not stained in Steps 2 or 3, aliquot 100 μ L of cell suspension to each tube. The cell number should be determined empirically but can range from 10^5 to 10^8 cells/tube.

For additional questions, please contact Technical Support at +1-888-810-6168 (US) or +43 1 796 4040 120 (Europe/International), or send us an email at Tech_Support@affymetrix.com

BrdU Staining Kit for Flow Cytometry

Research Use Only

- c) Gently resuspend the cells from Step 2f, Step 3d, or Step 4b, by pulse-vortexing once. This resuspension step is critical before the addition of the freshly prepared 1X BrdU Staining Buffer working solution.
- d) Add 1 mL of freshly prepared 1X BrdU Staining Buffer working solution and mix gently. Incubate for 15 minutes at room temperature in the dark. Incubations may go longer (up to 14 hours) but should be determined empirically for each cell type.
- e) Wash cells twice with Flow Cytometry Staining Buffer, as described in Step 1d.
- f) Add 100 μ L of the DNase I working solution that was prepared in Step 4a to each sample. Incubate for 1 hour at 37°C in the dark.
- g) Wash cells twice with Flow Cytometry Staining Buffer, as described in Step 1d.
- h) Add 5 μ L of Anti-BrdU fluorochrome-conjugated antibody per sample. Mix and incubate for 20-30 minutes at room temperature in the dark.
Note: Antibodies against intracellular antigens or surface antigens not stained in Step 3 may be added here. The antibodies used for surface staining at this step must recognize a fixed epitope. If an antibody only recognizes a native epitope or if this information is unknown, we recommend surface staining at Step 3.
- i) Wash cells twice with Flow Cytometry Staining Buffer, as described in Step 1d.

Step 5: Acquire data on a flow cytometer.

For additional questions, please contact Technical Support at +1-888-810-6168 (US) or +43 1 796 4040 120 (Europe/International), or send us an email at Tech_Support@affymetrix.com

Revised 08-09-16

© Affymetrix, Inc. All rights reserved

Tel: +1-888.999.1371 or +1-858.642.2058 • Fax: +1-858.642.2046 • ebioscience.affymetrix.com • info@ebioscience.com