

Revised: June 25, 2012

Product Information

Fixation Buffer

Catalog Number: 22015

Unit Size: 100 mL

Color and Form: Colorless solution

Storage and Handling: Store at room temperature. Do not freeze. Stable for at least 6 months from the date of receipt. Discard if cloudy.

Warning: Fixation Buffer contains formaldehyde, which is toxic by ingestion, inhalation, and contact with skin, and is a suspected carcinogen. Avoid contact with skin, eyes, and clothing. Dispose as toxic waste.

Product Description

Fixation Buffer is a ready-to-use formaldehyde-based fixative for immunofluorescence microscopy or flow cytometry.

Protocols

Fixation Buffer can be used in any standard cell fixation protocol. Protocols for Biotium's Fixation Buffer (22015), Permeabilization Buffer (22016), and Permeabilization and Blocking Buffer (22017) are provided below and may be optimized for specific applications.

Protocol for intracellular staining for flow cytometry

- 1. Pellet cells by centrifuging at 350 x g for 5 minutes. Wash cells twice in PBS. To wash cells, resuspend the cell pellet in PBS, centrifuge at 350 x g for 5 minutes, and gently pour off supernatant. Resuspend cells in PBS at a density of 10^7 cells/mL.
- 2. Aliquot 100 uL of cell suspension (10^6 cells) per tube into 12 x 75 mm polypropylene flow cytometry tubes.
- 3. Staining for surface antigens can be performed at this point:
 - a. Add the appropriate antibodies to cells in PBS.
 - b. Incubate for 15 minutes at room temperature in the dark.
 - c. Wash cells twice with 3 mL PBS as described in step 1.
 - d. Resuspend cells in 100 uL PBS.
- 4. Add 100 uL of Fixation Buffer to each tube and vortex gently to mix.
- 5. Incubate for 20 minutes at room temperature.
- Centrifuge for 5 minutes at 350 x g. Wash cells twice in PBS + 2% bovine serum or goat serum. To wash cells, resuspend cell pellet in 3 mL wash buffer, centrifuge for 5 minutes at 350 x g, and gently pour off supernatant.
- Add 100 uL of Permeabilization Buffer to each tube. Add primary antibodies to the permeabilization buffer at the antibody suppliers' recommended concentrations, and vortex gently to mix. A negative control omitting primary antibodies (or using isotype controls) is recommended to measure background.
- 8. Incubate at room temperature for 30 minutes. If staining with fluorescentlylabeled primary antibodies, incubate in the dark.
- 9. Wash cells twice with wash buffer (see step 6).
- If staining with fluorescently-labeled primary antibodies, add 1 mL wash buffer and analyze by flow cytometry. If staining with unconjugated primary antibodies and fluorescently-labeled secondary antibodies, proceed to next step.
- Resuspend the cells in the residual wash buffer remaining in the tube after step 9 (~100 uL). Add fluorescent secondary antibody conjugates at the suppliers' recommended concentrations and vortex gently to mix.

- 12. Incubate for 30 minutes at room temperature in the dark.
- 13. Wash cells twice with wash buffer (see step 6).
- 14. Resuspend cell pellet in 1 mL wash buffer and analyze by flow cytometry.

Immunofluorescence Protocol for Microscopy

Note: Because formaldehyde-based fixatives can compromise plasma membrane integrity, fixation at 4° C is recommended for staining of surface antigens in non-permeabilized cells.

1. Coverslip preparation for adherent cells

- 1.1 Culture cells on slide chambers or sterile glass coverslips (with poly-L-lysine coating if cells do not adhere well, see below). We recommend 18 x 18 mm square coverslips in 6-well plates or 4-well chamber slides.
- 1.2 Allow cells to adhere and culture or treat as desired.

2. Coverslip preparation for non-adherent cells

- 2.1 Coat coverslips with 0.01% poly-L-lysine solution for 10 minutes at room temperature.
- 2.2 Aspirate the poly-L-lysine solution and allow coverslips to dry completely.
- 2.3 Centrifuge cells in medium and resuspend in PBS. Transfer cells to coverslips.
- 2.4 Incubate for 30-60 minutes. Check for adherence under microscope.

3. Fixation and Staining

- 3.1 Rinse cells three times in PBS to remove culture medium.
- 3.2 Incubate cells in Fixation Buffer (catalog no. 22015), 15 minutes at room temperature. Alternatively, cells can be fixed in chilled (4°C) Fixation Buffer on ice for 30 minutes. Other fixatives also can be used.
- 3.3 Rinse twice with PBS to remove traces of fixative.
- 3.4 Incubate with 1X Permeabilization and Blocking Buffer, 30 minutes at room temperature.
- 3.5 Dilute primary antibody in 1X Permeabilization and Blocking Buffer to the concentration recommended by the antibody supplier. A negative control with primary antibody omitted is recommended to assess background. Overlay enough diluted antibody solution to completely cover cells. Parafilm® squares can be overlaid on coverslips to evenly spread a small volume (~100 uL) of antibody solution over the surface. Keep slips in a humidified chamber to avoid evaporation. Incubate 2 hours at room temperature, or overnight at 4°C.
- 3.6 Wash three times with PBS, 5 minutes each wash. 1X Permeabilization and Blocking Buffer can be used instead of PBS for a more stringent wash.
- 3.7 Dilute fluorescent secondary antibody in dilution buffer and incubate for 30 minutes to 2 hours at room temperature. IgG conjugates can be used at 1-10 μg/mL for most applications. Fluorescent nuclear stains or phalloidins can be included at this step.
- 3.8 Wash three times with PBS, 5 minutes each wash. 1X Permeabilization and Blocking Buffer can be used instead of PBS for a more stringent wash.
- 3.9 Mount coverslips using anti-fade mounting media, such as EverBrite™ Mounting Medium (see related products), or add enough mounting medium to wells or chambers to completely cover cells. Seal coverslip edges with clear nail polish or CoverGrip™ coverslip sealant (catalog no. 23005).
- 3.10 Store slides in the dark at 4°C.

Tips and Hints:

- No signal or weak fluorescence intensity may suggest the following: (a) insufficient antibody is present for detection, (b) intracellular target was not accessible, (c) excitation sources are not aligned, (d) target protein is not present or expressed at low levels, (e) fluorochrome has faded, and/or (f) primary and secondary antibodies are not compatible.
- High fluorescence background may suggest the following: (a) antibody concentration is too high, (b) excess antibody was not washed away efficiently, and/or (c) blocking was inadequate. Increase antibody dilution and washes.

Fixation Buffer

Related Products

Cat.#	Product Name	Unit Size
23006	Flow Cytometry Fixation/Permeabilization Kit	50 assays
22016	Permeabilization Buffer	100 mL
22017	Permeabilization and Blocking Buffer	100 mL
30069	AccuEasy™ Flow Cytometry Kit	1 kit
23001	EverBrite™ Mounting Medium	10 mL
23002	EverBrite™ Mounting Medium with DAPI	10 mL
23003	EverBrite™ Hardset Mounting Medium	10 mL
23004	EverBrite™ Hardset Mounting Medium + DAPI	10 mL
23005	CoverGrip™ Coverslip Sealant	15 mL
22010	10% Fish Gelatin Blocking Buffer	100 mL
22011	Fish Gelatin Powder	2 x 50 g
22014	30% Bovine Serum Albumin Solution	100 mL
22012	Dry Milk Powder	4 x 25 g
22002	Tween®-20	50 mL

Please visit www.biotium.com to view our full selection of products featuring bright and photostable fluorescent CF[™] dyes, including secondary antibodies and Mix-n-Stain[™] antibody labeling kits. Biotium also offers a variety of apoptosis and cell viability assays for flow cytometry analysis, including mitochondrial membrane potential dyes, fluorescent Annexin V conjugates, and NucView[™]488 Caspase-3 Substrate for live cells.

CF, Mix-n-Stain, and NucView are trademarks of Biotium. Tween is a registered trademark of Uniqema Americas LLC. Parafilm is a registered trademark of Bemis, Co.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.