

ReadiUse™ 4% formaldehyde fixation solution

Catalog number: 20010
Unit size: 50 mL

Component	Storage	Amount (Cat No. 20010)
ReadiUse™ 4% formaldehyde fixation solution	Refrigerated (2-8 °C), Minimize light exposure	1 bottle (50 mL)

OVERVIEW

All our ReadiUse™ reagents require minimal hands-on time. This 4% formaldehyde solution is a ready-to-apply format. It is widely used for fixing cells in fluorescence imaging applications.

AT A GLANCE

Important Note: Store at 4°C and avoid exposure to light.

SAMPLE EXPERIMENTAL PROTOCOL

Following is our recommended protocol which only provides a guideline. It should be modified according to your specific needs.

1. Perform formaldehyde fixation: Incubate the cells with 4% formaldehyde for 10–30 minutes at room temperature.

Note: Avoid any methanol containing fixatives since methanol can damage actin during the fixation process. The preferred fixative is methanol-free formaldehyde

2. Rinse the fixed cells 2–3 times in PBS.
3. The cells are ready for immunohistochemistry/immunofluorescence.

Note: Store the fixed cells in PBS at 4 °C.

EXAMPLE DATA ANALYSIS AND FIGURES

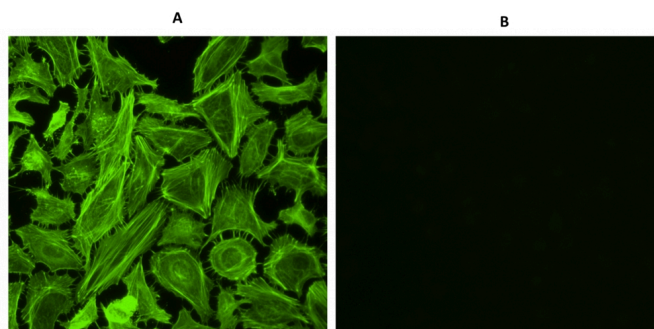


Figure 1. F-actin stain of CPA cells.

CPA cells were fixed with ReadiUse™ 4% formaldehyde fixation solution (Cat#20010) and then stained with Cell Navigator® F-Actin Labeling Kit (cat#22661). A: label with 1X iFluor® 488-phalloidin for 30 min only. B: Treat the cells with phalloidin for 10 min, then stain them with 1X iFluor® 488-phalloidin for 30 min.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.