Glowing products for science

## **Fixation Buffer**

Ready-to-use formaldehyde-based fixation buffer for immunofluorescence staining for microscopy or flow cytometry.



## **Product attributes**

Storage Conditions Store at room temperature

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

Fixation Buffer is a ready-to-use formaldehyde-based fixation buffer for immunofluorescence cell staining for microscopy or flow cytometry.

- Made with methanol-stabilized formaldehyde
- Fix cells for microscopy and flow cytometry applications
- 100 mL solution

The Fixation Buffer is also available in the Flow Cytometry Fixation/Permeabilization Kit (<u>23006</u>). This Fixation Buffer is prepared with methanol-stabilized formaldehyde. For a methanol-free formaldehyde fixative, see our <u>Paraformaldehyde</u>, <u>4% in PBS, Ready-to-Use Fixative</u>.

This datasheet was generated on November 13, 2025 at 10:25:32 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/fixation-buffer/">https://biotium.com/product/fixation-buffer/</a>