



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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

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 ([23006](#)). This Fixation Buffer is prepared with methanol-stabilized formaldehyde. For a methanol-free formaldehyde fixative, see our [Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative](#).

This datasheet was generated on June 9, 2026 at 08:43:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/fixation-buffer/>