



Glowing products for science

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

## Flow Cytometry Fixation/Permeabilization Kit

A kit containing fixation and permeabilization buffers used in staining intracellular antigens for flow cytometry.



### Product attributes

Storage Conditions
Store at room temperature

### Product Description

The Flow Cytometry Fixation/Permeabilization Kit contains optimally formulated buffers for fixation and permeabilization of cells for immunofluorescence staining of intracellular antigens for analysis by flow cytometry.

Note: the fixation buffer in this kit is prepared from methanol-stabilized formaldehyde. For methanol-free fixative, see our Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative ([22023](#)).

See our [Flow Cytometry Technology Page](#) to see all of our products for flow cytometry.

This datasheet was generated on January 1, 2026 at 06:02:17 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/flow-cytometry-fixationpermeabilization-kit/>