Glowing products for science

# **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

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

## **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

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 November 2, 2025 at 12:21:50 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/flow-cytometry-fixationpermeabilization-kit/