



Glowing products for science

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

Permeabilization and Blocking Buffer (5X)

A concentrated 5X buffer solution for one step permeabilization and blocking for immunofluorescence staining for fluorescence microscopy.



Product attributes

Storage Conditions

Store at -10 to -35 °C

Product Description

The Permeabilization and Blocking Buffer (5X) is a concentrated stock solution for one step cell permeabilization and blocking. It is designed for use during immunofluorescence staining for fluorescence microscopy.

- Permeabilizes fixed cells for immunofluorescence staining
- Blocks non-specific binding of antibodies
- Can be used to dilute antibodies for staining, and as a wash buffer
- Contains goat serum, not for use with anti-goat or anti-sheep antibodies
- 100 mL of 5X solution

Note: this buffer contains goat serum and therefore is not compatible with anti-goat and anti-sheep secondary antibodies. We recommend Permeabilization Buffer ([22016](#)) and 10% Fish Gelatin Blocking Agent ([22010](#)) for use with anti-goat and anti-sheep secondary antibodies.

This datasheet was generated on January 18, 2026 at 06:47:40 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/permeabilization-and-blocking-buffer-5x/>