

Permeabilization and Blocking Buffer (5X)

THE METERS OF THE PARTY OF THE

Storage Conditions

Call us: 800-304-5357

Product attributes

Store at -10 to -35 °C

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

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 November 22, 2025 at 11:08:34 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/permeabilization-and-blocking-buffer-5x/