Glowing products for science

Dry Milk Powder

A blocking agent commonly used to minimize background noise in immunodetection procedures such as western blotting and immunostaining.



Product attributes

Storage Conditions

Store at room temperature

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

Product Description

Dry Milk Powder can be dissolved to make blocking buffers to minimize non-specific antibody binding in immuno-detection procedures such as western blotting, immunofluorescence, and IHC.

- Nonfat milk
- Supplied as a dry powder
- Use for blocking during western blotting or immunostaining
- 4 x 25 g

Nonfat dry milk is commonly added to blocking buffers to minimize background noise in immunodetection procedures such as western blotting, immunofluorescence, and IHC. During western blotting, milk can be used to block non-specific binding to nitrocellulose or PVDF membranes. Milk can also be used for blocking and dilution of antibodies during immuno-staining of cells or tissues.

This item is only available for customers in the USA.

This datasheet was generated on November 9, 2025 at 06:20:04 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dry-milk-powder/