

# Product Information

## Non-fat Dry Milk Powder

**Catalog Number:** 22012

**Unit Size:** 4 X 25 g

**Color and Form:** White to light yellow powder

### Storage and Handling

Store at RT and protect from moisture.

### Product Description

Dry milk can be used as a blocking solution that can minimize background noise in immunodetection procedures such as Western Blotting and immunostaining. Milk blocking buffer and be incubated with membranes or for dilution of antibodies during immunostaining procedures.

### Protocol:

The protocol below is a general guideline. Some optimization may need to be required depending on the application.

### Reagents:

Reagents required, but not included:

- Tris Buffered Saline (TBS) or Phosphate Buffered Saline (PBS)
- Tween® 20 or other detergent of choice (optional)

#### 1. Preparation of 100 mL 1X blocking buffer:

Dissolve 1 - 5 g of dry milk in 100 mL PBS or TBS. Detergents such as Tween® 20 can be added to a final concentration of 0.1%.

Reagent	Volume
1X PBS for TBS	100 mL
Dry Milk	1 - 5 g
Tween 20 (optional)	100 µL

2. Incubate membrane in 1X Blocking Buffer for 30 to 60 minutes at room temperature with gentle agitation, or incubate overnight at 4°C with gentle agitation. Blocking for immunostaining can also be performed similarly on cells or tissues.
3. Continue Western blotting or staining procedure according to desired protocol. Procedures may need to be optimized and should be determined empirically.

### Related Products

Cat.#	Product Name	Unit Size
22010	Fish Gelatin Blocking Buffer, 10X	100 mL
22011	Gelatin, cold water fish	100 g
22013	BSA	50 g
22002	Tween 20	50 mL
23001	EverBrite™ antifade mounting medium	10 mL
23002	EverBrite™ antifade mounting medium with DAPI	10 mL

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.