

www.biotium.com

Revised: August 28, 2014

Product Information

10X Fish Gelatin Blocking Agent

Catalog Number: 22010

Unit Size: 100 mL

Storage and Handling

Store at 4°C. Product is stable for at least 6 months from date of receipt when stored as recommended. Warm to room temperature before use. 10X Fish Gelatin Blocking Agent solution will gel during storage at 4°C, it can be warmed to 37°C to liquify. Fish Gelatin Blocking Agent will not gel at 4°C after dilution to 1X in buffer.

Product Description

Gelatin purified from cold water fish skin is used as a blocking agent to reduce nonspecific antibody binding in immunodetection procedures such as Western Blotting and immunostaining. It is used as an alternative to BSA, milk, or mammalian serum because does not contain IgG that could cross-react with mammalian antibodies.

Protocol

General guidelines for using Fish Gelatin Blocking Buffer are provided below. Protocols may require optimization for specific applications. Blocking buffer containing 1X Fish Gelatin Blocking Agent can be used to replace blocking buffers containing 5% non-fat dry milk or 5% BSA.

Reagents required, but not provided

- Tris Buffered Saline (TBS) or Phosphate Buffered Saline (PBS)
- Tween® 20, Triton™ X-100, or other detergent of choice (optional)

Preparation of 1X Fish Gelatin Blocking Buffer

Warm 10X Fish Gelatin Blocking Agent to room temperature and gently swirl or invert bottle to mix well. Add 1 volume of 10X Fish Gelatin Blocking Agent to 9 volumes of TBS or PBS buffer for a final concentration of 1X Fish Gelatin Blocking Agent. Stir to mix thoroughly. Store blocking buffer at 4°C. Discard blocking buffer if it becomes cloudy.

Optional: A detergent such as Tween® 20, Triton™ X-100, or other detergent of choice may be added to blocking buffer at a final concentration of 0.1%.

Optional: A preservative such as sodium azide (0.02% final concentration) may be added to blocking buffer for long term storage (note: sodium azide is not compatible with horseradish peroxidase (HRP) based assays).

Blocking and staining

Incubate Western membranes in 1X Blocking Buffer for 30 to 60 minutes at room temperature with gentle agitation. Blocking for immunostaining can also be performed similarly on cells or tissues. 1X Blocking Buffer can be used as an antibody diluent for antibody incubation steps, and can be used during wash steps to reduce background.

Note: After dilution to 1X in buffer, Fish Gelatin Blocking Agent will not gel during incubation at 4°C.

Related Products

Catalog No.	Product Name
22011	Fish gelatin powder
22013	Bovine serum albumin, Fraction V
22014	Bovine serum albumin, 30% solution
22012	Non-fat dry milk
22002	Tween® 20
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
23005	CoverGrip™ Coverslip Sealant
22005	Mini Super ^{н⊤} Pap Pen 2.5 mm tip, ~400 uses
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent and near-infrared CF™dye antibody conjugates and antibody labeling kits, HRP, R-PE, and APC antibody conjugates and antibody labeling kits, plus apoptosis reagents, qPCR reagents, and other tools for cell biology research.

Tween is a registered trademark of Uniqema Americas LLC. Triton is a trademark of Dow Chemical Company.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.