

Cat. # T7131A

For Research Use

TAKARA

**Western BLoT Blocking Buffer
(Fish Gelatin)**

Product Manual

v201403Da_201802

Table of Contents

I. Description	3
II. Components	3
III. Storage	3
IV. Precautions.....	3
V. Protocol.....	3
VI. Related Products	4

I. Description

The Western BLoT Blocking Buffer (Fish Gelatin) is a blocking buffer that contains gelatin extracted from fish in TBS-T buffer. Since the blocking reagent in this product is of fish origin, it is less likely to cross react with antibodies of mammalian origin than conventional blocking reagents of mammalian origin, such as dry milk and BSA. With the low background obtained with this product, it is possible to clearly detect antigens.*

This Blocking Buffer is ready-to-use without dilution for blocking Western blot membranes. The Buffer can also be used for dilution of primary antibodies and secondary antibodies labeled with AP or HRP.

- * : There is no common blocking reagent for Western blotting that can be used for all experiments. Preliminary testing is necessary to determine the optimal blocking reagent. Depending on the specific antigen, blocking with gelatin may decrease detection sensitivity because the antigen may be masked by gelatin. In this case, a better result may be obtained with another blocking reagent, such as BSA or casein.

II. Components

Western BLoT Blocking Buffer (Fish Gelatin) 1 L

III. Storage 4°C

IV. Precautions

The following are precautions for using this product. Read before use.

1. After using the Blocking Buffer, close the lid immediately to prevent contamination.
2. Bring the Blocking Buffer to room temperature before use.
3. The optimal blocking conditions depend on the experimental system. Determine the optimal blocking time at room temperature.

V. Protocol

1. After protein is transferred to the Western blot membrane, completely cover the membrane with an appropriate volume of Western BLoT Blocking Buffer (Fish Gelatin). Shake gently for 10 - 60 minutes at room temperature. The membrane can also be blocked overnight at 4°C. Use approximately 0.2 ml of the buffer per cm² of membrane.
2. After blocking, perform detection with appropriate antibodies. This product can also be used for dilution of primary and secondary antibodies.

VI. Related Products

Western BLoT Chemiluminescence HRP Substrate (Cat. #T7101A/B)
Western BLoT Quant HRP Substrate (Cat. #T7102A/B)
Western BLoT Hyper HRP Substrate (Cat. #T7103A/B)
Western BLoT Ultra Sensitive HRP Substrate (Cat. #T7104A/B)
Western BLoT Immuno Booster (Cat. #T7111A)
Western BLoT Rapid Detect (Cat. #T7121A)
Tris-Glycine-SDS Buffer (TG-SDS) Powder, pH8.3 (Cat. #T9101)
Tris-Glycine Buffer (TG) Powder, pH8.3 (Cat. #T9102)
CLEARLY Protein Ladder (Unstained) (Cat. #3453A/B)
CLEARLY Stained Protein Ladder (Cat. #3454A/B)
Phosphate Buffered Saline with Tween20 (PBS-T) Tablets, pH7.4 (Cat. #T9183)
Tris Buffered Saline with Tween20 (TBS-T) Tablets, pH7.6 (Cat. #T9142)
Western BLoT Stripping Buffer (Cat. #T7135A)

NOTE: This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc.

Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from TAKARA BIO INC.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at www.takara-bio.com.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.
