WesternBreeze® Chemiluminescent Western Blot Immunodetection Kit



Package Contents

Catalog Numbers WB7104 Anti-Mouse WB7106 Anti-Rabbit

WB7108 Anti-Goat





- Store all solutions at 4°C.
- Once diluted, use the solutions the same day.
- All solutions are proprietary formulations and contain chlorobutanol as a preservative.



Required Materials





Timing

Preparation and immunodetection: ~3 hours



Selection Guide

Western Blotting Kits

Go online to view related products.



Product Description

- WesternBreeze® Chemiluminescent Kits detect proteins that have been immobilized on membranes (nitrocellulose or PVDF) following western transfer or bound directly from solution (dot blots).
- Detection is accomplished with a ready-to-use CDP-Star® chemiluminescent substrate for alkaline phosphatase.
- Protein bands can be captured either by X-ray film or a CDP-Star® compatible imaging system.



Important Guidelines

- Avoid touching the working surface of the membrane, even with gloves.
- Use ultra-filtered water, free from alkaline phosphatase activity. Do not allow the membrane to dry out after adding the chemiluminescent substrate.
- Use a rotary shaker platform rotating at 1 revolution/ second for all washing, blocking, and incubating steps.
- Add solutions to the trays slowly, at the membrane edge, to avoid bubbles forming under the membrane. Decant from the same corner of the dish to ensure complete removal of previous solutions.



Online Resources

Visit our product pages for additional information and protocols. For support, visit www.lifetechnologies.com/support.





Protocol Outline

This process involves the following phases:

- A. Prepare membrane and solutions.
- B. Incubate membrane with Primary and then Secondary Antibodies.
- C. Add and incubate substrates.
- D. Complete detection.

Western Blotting Protocol

- **1** See page 2 for quantities and guidelines to prepare the antibodies and solutions.
- 1 See page 3 to view a typical immunodetection procedure for use with small membranes (60 cm²).

Preparing the Membrane

For Westerns or Western dot blots, wash membrane twice for 5 minutes in 20 mL of water to remove gel, transfer buffer, and weakly bound proteins. Then, proceed with immunodetection.

For water washed and dried membranes:

- NC Membranes: Proceed with immunodetection.
- PVDF membranes: Re-wet membrane in methanol followed by two 20-mL water washes for 5 minutes.

Scaling up for Large Membranes

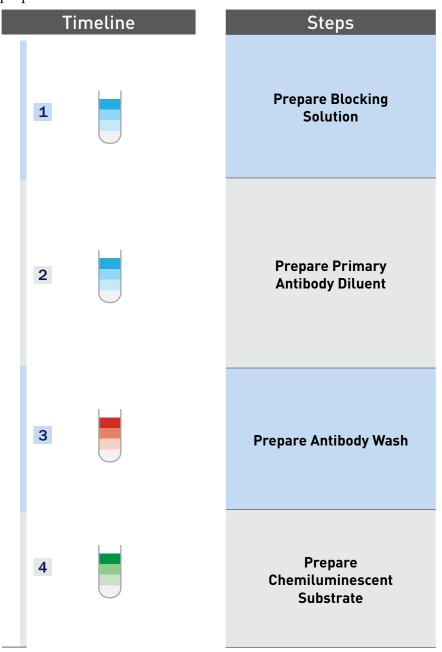
To blot standard size gels (\sim 200 cm²), scale up the required solution volumes \times 3.3. Use a tray that closely matches the dimensions of the membrane for the most efficient use of the solutions.



For Research Use Only. Not for use in diagnostic procedures.

Preparing Solutions for Nitrocellulose and PVDF Membranes

The Blocking Solution is used for blocking and as a Primary Antibody Diluent for NC membranes. For western blots from Novex $^{\circ}$ or other mini gels, prepare solutions as described in table below for $\sim 60 \text{ cm}^2$ membrane.



Procedure Details					
For NC Membran	е	For PVDF Membrane			
Ultra filtered Water	14 mL	Ultra filtered Water	5 mL		
Blocker/Diluent (Part A)	4 mL	Blocker/Diluent (Part A)	2 mL		
Blocker/Diluent (Part B)	2 mL	Blocker/Diluent (Part B)	3 mL		
Total Volume	20 mL	Total Volume	10 mL		
For NC Membran	е	For PVDF Membrane			
Dilute your Primary Antibody according to the manufacturers recommendations into 10 mL of NC Blocking Solution (see above). Typically, commercial Primary Antibody preparations are diluted 1:1000 to 1:5000 to a concentration of about 1 to 0.2 µg/mL.		Ultra filtered Water	7 mL		
		Blocker/Diluent (Part A)	2 mL		
		Blocker/Diluent (Part B)	1 mL		
		Total Volume	10 mL		
		Dilute your Primary Antibody into this diluent according to manufacturer's recommendations.			
For NC Membran	е	For PVDF Membrane			
Ultra-filtered Water	150 mL	Ultra-filtered Water	150 mL		
Antibody Wash Solution (16X)	10 mL	Antibody Wash Solution (16X)	10 mL		
Total Volume	160 mL	Total Volume	160 mL		
For NC Membrane		For PVDF Membrane			
Chemiluminescent Substrate	2.375 mL	Use the CDP-Star directly from the bottle. DO NOT add the Chemilumi-			
Chemiluminescent 0.125 mL Substrate Enhancer		nescent Substrate enhancer (Nitro-Block-II enhancer).			
Total Volume	2.5 mL	Total Volume	2.5 mL		

Immunodetecting with the Chemiluminescent Western Blot Kit

The table below describes the procedure for immunodetection using small membranes (60 cm²).

	T	imeline	Steps	Procedure Details
	1		Block membrane	Use 10 mL of blocking solution in the dish provided. Incubate 30 minutes on a rotary shaker (1 revolution/second). Decant the Blocking Solution.
	2		Wash membrane with water	Rinse the membrane with 20 mL of water for 5 minutes, then decant. Repeat once.
	3		Incubate with Primary Antibody	Incubate the membrane with 10 mL of Primary Antibody Solution for 1 hour, then decant.
	4		Wash membrane with Antibody Wash	Wash membrane for 5 minutes with 20 mL of prepared Antibody Wash, then decant. Repeat 3 times.
	5		Incubate with Secondary Antibody	Incubate the membrane in 10 mL of Secondary Antibody Solution for 30 minutes, then decant.
Day 1	6		Wash membrane with Antibody Wash	Wash the membrane for 5 minutes with 20 mL of Antibody Wash, then decant. Repeat 3 times.
	7		Wash membrane with water	Rinse the membrane with 20 mL of water for 2 minutes, then decant. Repeat twice.
	8		Add substrate	Place the membrane on sheet of transparency plastic. Do not allow membrane to dry out. Evenly apply 2.5 mL of Chemiluminescent Substrate Solution to the membrane without touching the surface (remember to add Chemiluminescent Substrate Enhancer for NC membranes).
	9	5	Incubate with substrate	Allow reaction to develop for 5 minutes at room temperature.
	10		Detect	Blot excess Chemiluminescent Substrate from membrane using filter paper in the kit. Do not allow membrane to dry out. Cover membrane with another piece of transparency plastic. Expose x-ray film. for one second to several minutes, or use another imaging device.