

Protocol Using TidyBlot Western Blot Detection Reagent

WB

For use with TidyBlot HRP Conjugated Western Blot Detection Reagent. TidyBlot Reagent is very versatile because it is compatible with a variety of both monoclonal and polyclonal primary antibodies. [View TidyBlot's compatibility chart.](#) TidyBlot preferentially binds to the native antibody used during the western blotting procedure rather than the SDS-denatured/reduced form present in lysates/immunoprecipitates. Best results are therefore achieved by denaturing the immunoprecipitate/lysate fully. If your target protein is phosphorylated, we recommend using our [protocol for detection of phosphorylated proteins](#) by western blotting. Please note that a certain level of technical skill and immunological knowledge is required for the successful design and implementation of these techniques - these are guidelines only and may need to be adjusted for particular applications.

Reagents

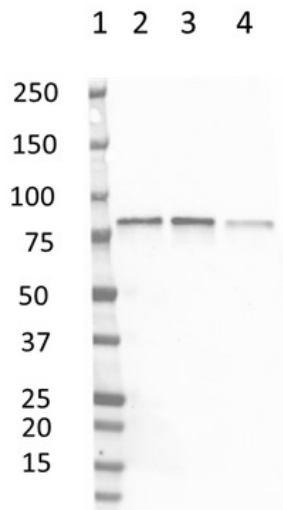
- **Blocking Buffer**
Block Ace (Cat. #BUF029) dissolved in ddH₂O (100 ml of ddH₂O per sachet of Block Ace)
- **TGS Running Buffer**
100 ml 10x TGS Running Buffer (#161-0772) diluted in 900 ml deionized water (DI H₂O)
- **Washing Buffer:**
Block Ace (#BUF029) dissolved in PBS (1 L of PBS per sachet of Block Ace) +0.1% Tween 20
- **Ponceau S**
Acetic acid, 5 ml
Distilled water, 95 ml
Ponceau S (Sigma, #P3504), 0.1 g
*Note: Ponceau S is light sensitive.

Method

1. Prepare samples of your protein of interest in sample buffer such as 2x Laemmli Sample Buffer (#1610737) with reducing agent, for example Dithiothreitol (DTT) (#1610610). Boil the samples for 5 min at 95 °C and allow to cool to room temperature (RT). Please note that this step is vital as complete reduction and denaturation of proteins in your samples is required to prevent TidyBlot Reagent from detecting antibodies present in your samples.
2. Place an SDS page gel into a suitable gel tank. The size and type of gel selected should be optimized for the system you are using and the size of your target protein. In general, gels with a higher percentage of acrylamide are more suitable for targets of a lower molecular weight.
3. Fill the inner chamber and outer chamber of the gel tank with running buffer according to manufacturer's instructions.
4. Load equal amounts of prepared lysate into the wells of an SDS page gel. Remember to load a suitable molecular weight marker such as Bio-Rad's Precision Plus All Blue Standards (#161-0373) alongside your samples to determine the molecular weight of any proteins you visualize at the end of the blotting process.
5. Perform gel electrophoresis using a suitable protocol. For Bio-Rad's Criterion Gel System, we recommend running the gel at 300 V for approximately 25 min at RT.
6. Following SDS-PAGE, transfer proteins onto blotting membrane using a suitable transfer system, such as Bio-Rad's Trans-Blot Turbo Transfer System, according to the manufacturer's instructions.
7. Once transfer is complete, check that proteins have transferred from the gel to the membrane efficiently by staining the membrane with a total protein stain. Our recommendation is to stain the membrane post transfer with Ponceau S for 1 min, then completely destain the blot by washing with DI H₂O with agitation.
8. Place the membrane into blocking solution and incubate for 30 min at RT with agitation.
9. Remove the blocking buffer and wash the blot in wash buffer for 5 min with agitation.
10. Remove the wash buffer and add the primary antibody diluted in the wash buffer. Incubate overnight at 4°C with agitation. Please consult the product specific datasheet for the primary antibody you are using for recommended dilutions, bearing in mind that titration experiments may need to be performed in order to find the optimal primary antibody dilution for your experimental set-up.
11. Remove the primary antibody. Wash the blot in wash buffer 3x for 10 min with agitation.

12. Remove wash buffer and add TidyBlot Reagent diluted in wash buffer. We recommend diluting it between 1/40 and 1/400 in wash buffer. However, titration experiments should be performed to optimize the dilution of TidyBlot for the system you are using and your target protein. Incubate for 1 hr at RT with agitation.
13. Wash the membrane 4x for 5 min in wash buffer with agitation.
14. Wash the membrane 2x for 5 min in PBS.
15. Add an appropriate enzyme substrate solution to the membrane, such as Bio-Rad's Clarity ECL (#170-5061), and incubate as recommended by the manufacturer.
16. Develop the blot using an appropriate developing system for example Bio-Rad's ChemiDoc MP Imager according to manufacturer's instructions.

Appropriate controls should always be carried out. It may be useful to include a sample in which no primary antibody is used at all, in order to determine any nonspecific binding of the secondary reagent to the target tissue. Please contact [Bio-Rad's Technical Services Department](#) to learn about recommended secondary reagents for specific applications.



Western blot image using TidyBlot HRP conjugated Western Blot Detection Reagent. Western blot analysis of Protein Kinase C alpha (#MCA1572). A 25 µg amount of A549, Hepg2, and K562 cell lysates were loaded into lanes 2, 3, and 4 respectively. These cell lysates were run under reducing conditions and transferred onto a nitrocellulose membrane. Bio-Rad Precision Plus All Blue Standards (molecular weight marker) were run in lane 1. Mouse Anti-Protein Kinase C alpha (#MCA1572) was used as the primary antibody at a dilution of 1/1,000 in lanes 1-4. Goat Anti-Mouse IgG conjugated to HRP (#STAR207P) was used as the secondary reagent at 1/10,000. Membranes were visualized on the ChemiDoc MP Imaging System.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad-antibodies.com