

Avoiding Housekeeping Protein Detection Saturation

Protocol

Bulletin 6359

Reliable western blot data require detection of the target and loading control proteins in the linear dynamic range. Many published western blot images show saturated protein bands, especially for the housekeeping proteins (HKP) including β -actin, β -tubulin, and GAPDH, indicating that they were not measured in the linear dynamic range. Quantitative analysis based on this type of data is not reliable. This protocol describes common ways of avoiding signal saturation for the detection of HKP loading controls.

Method 1: Decrease sample load for HKP detection

Utilizes two gels to ensure quantitative detection of both target and HKP loading control proteins. One gel is used for target protein detection using a high sample load. A second gel is used for HKP detection using a very small amount of cell lysate to ensure the HKP detection is in the linear dynamic range.

Protocol:

1. Prepare two identical gels.
2. Load 20 μ g cell lysate for each sample in gel 1 for target protein detection.
Note: Many protocols use 10–50 μ g of cell lysate for the detection of target proteins. The appropriate amount of protein load should be determined before an actual experiment is performed. Please refer to *Determining the Appropriate Sample Load for Western Blots*, bulletin 6362, for details.
3. Load 1 μ g of cell lysate for each sample in gel 2 for HKP detection (such as β -actin and β -tubulin).
Note: The samples should be loaded in the same order as in gel 1. The appropriate amount of protein load should be determined before an actual experiment is performed. Please refer to *Determining the Appropriate Sample Load for Western Blots*, bulletin 6362, for details. A reasonable starting point is 1 μ g.
4. To ensure similar transfer conditions for both gels, use the same tank for a wet transfer, or the same cassette if using a fast transfer system.
5. Following the antibody manufacturer's instructions, apply the primary and secondary antibodies for detection of the target and loading control proteins on blots 1 and 2.
6. Develop the blots with a substrate, preferably one with a long signal duration such as Clarity™ western ECL substrate.
7. Use a CCD camera-based imager to capture the chemiluminescent signals on both blots. It is not recommended to use film.
8. The chemiluminescent image of blot 1 reveals the change in target protein expression level in different samples. The image of blot 2 reveals the load difference between the samples and can be used to normalize the target protein signals ensuring that the change observed in blot 1 is a true response to a biological event.

Note:

- Procedure details are omitted for steps such as protein sample prep, gel load, gel electrophoresis, transfer, or antibody incubation. For details, please refer to the *General Protocol for Western Blotting*, bulletin 6376
- Some journal reviewers may critique the fact that the loading control is from a different blot in a western blot experiment (Neill, U 2009)
- Once the experimental setup and conditions are established for the assay, do not change the sample load, transfer method, transfer time, antibody dilution, antibody incubation time, or temperature in subsequent experiments, as these factors may significantly change the detection signals

Method 2: Titration of antibody concentration

Utilizes only one gel for both the target and loading control protein detection. If the target is a low abundance protein and difficult to detect, the sample load is frequently increased in order to detect the target protein. It is common to titrate down the HKP antibody concentration to avoid saturation and achieve quantitative detection.

Protocol:

1. Select a typical sample in which an average amount of target and loading control proteins are present.
2. Load 20 μ g of the sample in each lane of a 15-well gel. Separate every two sample lanes with a prestained protein standard. Utilize two of the wells to load 1 μ g of the sample as reference.
Note: Typically, 10–50 μ g of cell lysate are used to detection target proteins. The appropriate amount of protein load should be determined before a real experiment is performed. Please refer to *Determining the Appropriate Sample Load for Western Blots*, bulletin 6362, for details.
3. Select a wet transfer or fast transfer method to transfer the gel. Following the transfer, cut the membrane along the prestained standard lanes to generate five blots.

BIO-RAD

4. Prepare the primary antibody against the housekeeping protein (e.g. β -actin) at different dilutions:
 - a. Manufacturer recommended dilution (e.g. 1:1,000)
 - b. 5x more dilute than recommended (e.g. 1:5,000)
 - c. 10x more dilute than recommended (e.g. 1:10,000)
 - d. 20x more dilute than recommended (e.g. 1:20,000)
5. Apply the four different dilutions of HKP primary antibody to the four blots loaded with 20 μ g of cell lysate. Treat the reference blot loaded with 1 μ g of cell lysate with the primary antibody using the manufacturer's recommended concentration.
6. Develop the blots using a substrate with a very long signal duration, for example, Clarity western ECL substrate.
7. Use a CCD camera-based imager to capture the chemiluminescent signals. It is not recommended to use film.
8. Based on the results from the blots, select the antibody concentration that most closely matches the data for the 1 μ g blot.

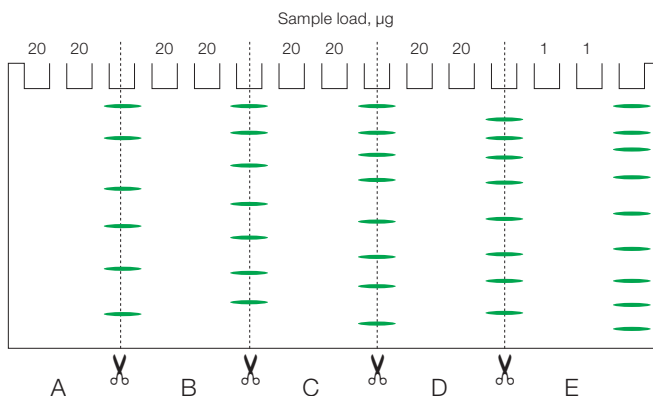
Note:

- Procedure details are omitted for steps such as protein sample prep, gel load, gel electrophoresis, transfer, or antibody incubation. For details, please refer to the *General Protocol for Western Blotting*, bulletin 6376
- Once the experimental setup and conditions are established for the assay, do not change the sample load, transfer method, transfer time, antibody dilution, antibody incubation time, or temperature in subsequent experiments, as these factors may significantly change the HKP signals

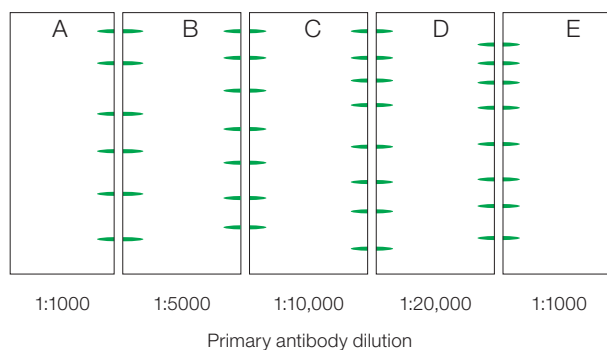
Reference

Neill U (2009). All data are not created equal. *J Clin Invest* 119, 424.

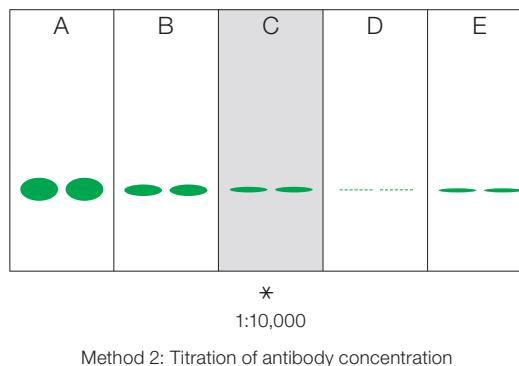
Gel Electrophoresis and Transfer



Antibody Incubation



Imaging and Analysis



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000