

Determining the Appropriate Sample Load When Performing a Stain-Free Western Blot

Protocol

Western Blotting

Bulletin 6363

Abstract

Traditional western blot workflows require an appropriate sample load to detect both target proteins and loading control proteins. The discrepancy between these quantities often leads to oversaturated protein bands, forcing researchers to sometimes run duplicate gels or titrate down the antibody concentration in order to get quantitative results.

Stain-Free western blotting requires only a sample load appropriate for target protein detection. No reprobing of the blot for loading control proteins is needed. More importantly, the Stain-Free blots exhibit greater linearity in the load range of 10–50 μg of cell lysate.

This protocol describes an assay development step to determine the right load for the detection of target proteins before the actual western blot experiment is carried out.

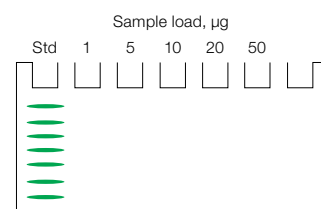
Protocol

1. Select a typical sample in which the target and loading control proteins are presented at an average level.
2. Load 1, 5, 10, 20, and 50 μg of protein sample to each lane.
3. After transfer, apply the target protein primary antibody to the blot.
4. Add substrate to develop the chemiluminescent signal and capture the signals using a CCD camera-based imager.
5. Use software to read the intensity of the target protein in each lane and plot those intensities against the protein load. Determine the linear dynamic range under the specific experimental conditions. Select the protein load that gives the quantitative reading in the linear dynamic range for the real experiment.

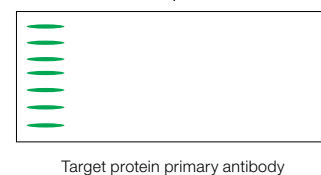
Note

- Procedure details are omitted for steps such as protein sample preparation, gel loading, gel electrophoresis, transfer, antibody incubation, etc. 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 detection signals

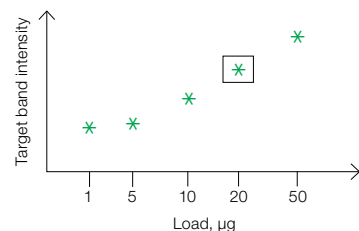
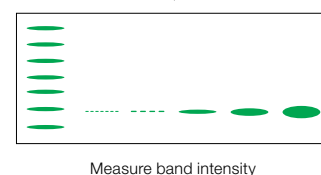
Gel Electrophoresis and Transfer



Antibody Incubation



Imaging and Data Analysis



BIO-RAD

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



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300
Israel 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23
Mexico 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23
Portugal 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23
Spain 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311
United Arab Emirates 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

