

Protocol

Determining the Linear Range for Quantitative Western Blot Detection



Published March 2017. Revised September 2019. The most recent version of this document is posted at [licor.com/bio/support](https://www.licor.com/bio/support).

Visit us on protocols.io! Explore an interactive version of this protocol at bit.ly/LinearRange-Overview.

Table of Contents

	Page
I. Introduction	2
II. Linear Range, Saturation, and Proportional Signals	4
III. Key Factors	5
IV. Required Reagents	5
V. Protocol: Determining the Linear Range for a Target Protein and Revert™ 700 Total Protein Stain	6
Total Protein and Target Quantification	7
Total Protein Quantification	7
Target Protein Quantification	8
Estimate Combined Linear Range for Target and Total Protein	9
VI. Protocol: Determining the Linear Range for a Target Protein and HKP	11
Target Protein and HKP Quantification	12
HKP Quantification (700 nm)	12
Target Protein Quantification (800 nm)	13
Estimate Combined Linear Range for Target and HKP	15
VII. Further Reading	17
References	17

I. Introduction

In quantitative Western blotting (QWB), normalization mathematically corrects for unavoidable sample-to-sample and lane-to-lane variation by comparing the target protein to an internal loading control. The internal loading control is used as an indicator of sample protein loading, to correct for loading variation and confirm that changes observed in target protein bands represent actual differences between samples.

QWB analysis is accurate only if the target protein and internal loading control can both be detected within the same linear range – a range that must be determined experimentally for each target and loading control. The combined linear range is then used to determine how much sample should be loaded to produce a linear signal response for both the target protein and the internal loading control (Figure 1).

More Info: Empiria Studio® Software provides a dedicated workflow for this process. Go to licor.com/empiria to learn more.

II. Linear Range, Saturation, and Proportional Signals

In QWB analysis, the **linear range of detection** is the range of sample loading that produces a linear relationship between the amount of target on the membrane and the band intensity recorded by the detector (Figure 1). Outside this range, signal intensity is not dependent on sample loading and does not accurately reflect the amount of target.

Within the linear range of detection, **band intensity should be proportional to the amount of target**. A change in target abundance should produce an equivalent change in signal response. For example, a two-fold increase in sample loading is expected to produce a two-fold increase in band intensity. At the upper and lower ends of the linear range, this proportional relationship is lost. Band intensity no longer reflects the abundance of target, and quantification is not possible.

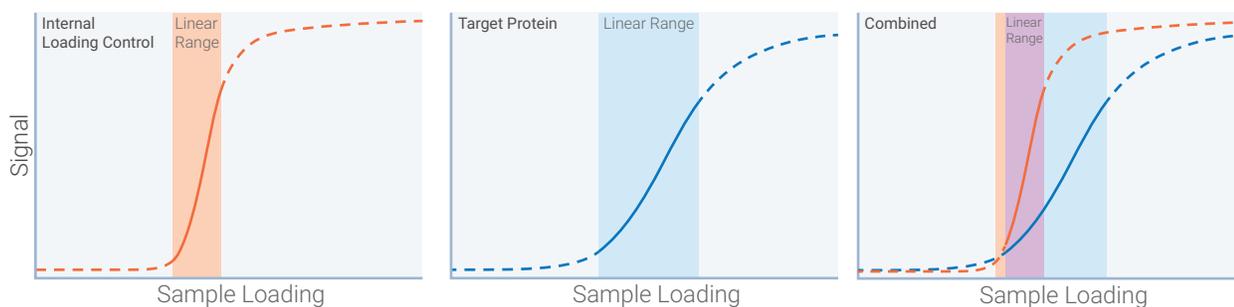


Figure 1. In the linear range of detection, the target and internal loading control both display a linear relationship between sample loading and band intensity. Determine the linear range separately for the target and internal loading control. Combine these ranges to identify a level of sample loading that will produce a linear response for both. Dotted lines indicate loss of proportionality on the upper and lower ends of the linear range.

Saturation occurs when increasing amounts of target fail to produce the expected increase in band intensity. Saturated bands yield similar intensity values, regardless of the actual amount of target present, and cannot be accurately quantified. Intensity of strong bands is underestimated, interfering with comparison of relative protein levels across the blot. Error introduced by saturation may alter data analysis and interpretation.

- **Membrane saturation** is the result of sample overloading. It frequently interferes with accurate detection of abundant proteins, including HKP loading controls. When samples are overloaded, abundant proteins can bind in layers on the membrane surface that limit antibody access during detection. Highly abundant proteins may also exceed the local binding capacity of the membrane and be washed away.
- **Signal saturation** occurs when the signal intensity of a band is too bright for the detection system to record. Increasing amounts of target do not produce a proportional increase in the recorded signal.

This protocol explains how to use serial dilutions of sample protein to determine the linear ranges of detection for a target and internal loading control, and choose an appropriate amount of sample to load for QWB analysis.

Note: This protocol is intended for use with near-infrared fluorescent Western blots.

III. Key Factors

Linear range of sample loading. This range must be determined individually for the target protein and internal loading control. Compare the linear ranges to determine the amount of sample you should load to produce a linear response for both target and control.

Housekeeping proteins. With two proteins of very different abundance, such as a target protein and highly expressed HKP loading control, it is absolutely critical to detect both proteins within the linear range. If the HKP is not detected in the linear range of sample loading, these strong bands will be underestimated and normalization will not be accurate.

Quantification of faint bands. Faint bands that are near the lower limit of detection generally have much larger coefficients of variation. They are difficult to discriminate from membrane background and may not be statistically significant. The high variability of faint bands introduces error that may affect data analysis and interpretation.

Antibody validation. Two-color Western blot detection requires careful selection of primary and secondary antibodies to prevent cross-reactivity. Always perform single-color control blots first to verify antibody specificity, and to identify possible interference from background bands.

IV. Required Reagents

1. **Treated and untreated samples**

Samples should represent the specific range of treatments or conditions you will use in your experiment (drug treatment, time course, dose response, etc.). Protein concentration must be determined for all samples.

2. **Revert™ 700 Total Protein Stain Kit (licor.com/revertkit)**

Revert 700 Total Protein Stain is used to assess sample protein loading in each lane as an internal loading control. After transfer and prior to immunodetection, the membrane is treated with this near-infrared fluorescent protein stain and imaged at 700 nm. Membrane staining can verify that sample protein was uniformly loaded across the gel, and assess the quality and consistency of protein transfer.

3. **Electrophoresis reagents**

4. **Transfer reagents**

5. **Western blot detection reagents (near-infrared fluorescence)**

Perform near-infrared Western blot detection according to the *Near-Infrared Western Blot Detection Protocol* (licor.com/NIRWesternProtocol; LI-COR).

V. Protocol: Determining the Linear Range for a Target Protein and Revert™ 700 Total Protein Stain

More Info: Empiria Studio® Software provides a dedicated workflow for this process. Go to licor.com/empiria to learn more.

Follow the instructions in this section if total protein staining of the membrane will be used as the internal loading control for QWB normalization. For normalization with a housekeeping protein, see Section VI.

1. Perform a Western blot with a serial dilution of sample (an 8-12 point, two-fold serial dilution is a good place to start).

Note: Make sure the dilution series includes sample loading amounts above and below the amount of sample you expect to load in each lane.

2. Prepare samples to be loaded on the gel with sample loading buffer.
3. Denature sample by heating at 95 °C for 3 min (or 70 °C for 10 min).
4. Load a uniform amount of sample protein in each lane.
5. Separate sample proteins by SDS-PAGE.
6. Transfer proteins to blotting membrane.
7. After transfer is complete, stain membrane with Revert™ 700 Total Protein Stain using the Single-Color Western Blot protocol:

More Info: For more information about the Revert protocol, see licor.com/revert.

- a. Add methanol to the stain reagents as indicated on each bottle.
- b. Rinse membrane in water, and incubate in 5 mL of Revert 700 Total Protein Stain solution for 5 minutes, with gentle shaking.
- c. Decant Total Protein Stain solution thoroughly. Rinse membrane twice with Wash solution.

Revert Wash solution (P/N 926-11012): 6.7% (v/v) Glacial Acetic Acid, 30% (v/v) Methanol, in water.

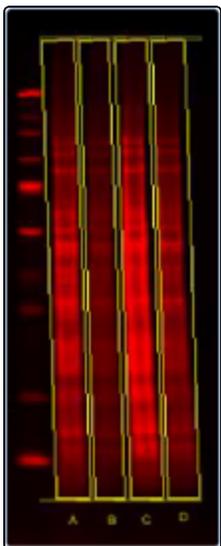
8. Image the membrane in the 700 nm channel with an Odyssey[®] imaging system (Classic, CLx, or Fc).
Adjust settings so that no saturation appears in the bands to be quantified.
9. Rinse membrane briefly with water, and proceed immediately to blocking and immunodetection steps.
10. Perform immunoblotting using IRDye[®] 800CW conjugated secondary antibody to detect the target protein.
11. Image the membrane in the 800 nm channel with an Odyssey imaging system.
Adjust settings so that no saturation appears in the bands to be quantified.

Total Protein and Target Quantification

Use Image Studio™ Software (licor.com/islite) to quantify the fluorescent signals from Revert staining (700 nm) and your target protein (800 nm).

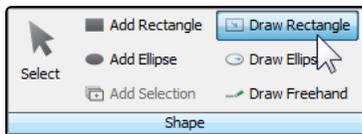
Total Protein Quantification

Use the Draw Rectangle tool in Image Studio™ Software to quantify the Total Protein signal in each lane.



1. Draw a shape around the first lane on the Total Protein Stain image (700 nm channel image).

a. In the Shape group on the Analysis tab, click Draw Rectangle.



b. Draw a rectangle around the first lane.
If lanes are skewed, rotate the shape by clicking Rotate in the Edit group.

More Info: For help choosing the right background subtraction method, see licor.com/BgSubtractHelp.

2. Add the shape from the first lane to each remaining lane.

- a. With the first shape selected, click Add Selection.
- b. Click the next lane to copy the shape.
If necessary, adjust the shape borders to ensure that each shape encloses the signal from one, and only one lane.

3. Export total protein quantification data.

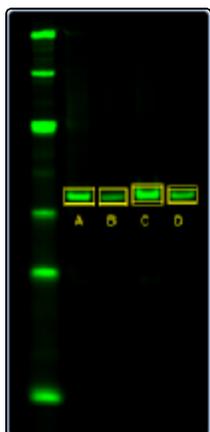
a. Click Shapes to open the Shapes data table.



b. Select shape data, then copy and paste data into a spreadsheet.
All data fields will be exported, but "Signal" is the field of interest for analysis.

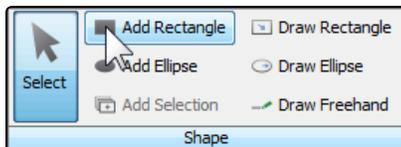
Target Protein Quantification

Use the Add Rectangle tool in Image Studio Software to quantify the target protein band in each lane (800 nm channel).



1. Click to select your Western blot 800 channel image, then add shapes to bands.

- a. In the Shape group on the Analysis tab, click Add Rectangle.



- b. Click each band to be analyzed, and an appropriately sized shape will be added around the band.

More Info: For help choosing the right background subtraction method, see licor.com/BgSubtractHelp.

2. Export quantification data.

- a. Click Shapes to open the Shapes data table.



- b. Select shape data, then copy and paste data into a spreadsheet.

Note: All data fields will be exported, but "Signal" is the field of interest for analysis.

Estimate Combined Linear Range for Target and Total Protein

1. Plot signal intensity vs. sample loading for both the target and Revert™ 700 Total Protein Stain (as shown in Figure 2).
2. Estimate the boundaries of the linear range of sample loading for the target and the total protein stain (shaded regions, Figure 2).
 - a. Estimate the low end of the linear range by selecting a level of sample loading that is above the level of membrane background (black dashed lines, Figure 2).
 - b. Estimate the upper end of the linear range by selecting a level of sample loading that is below the region where band intensity begins to plateau due to saturation (red dashed lines, Figure 2).

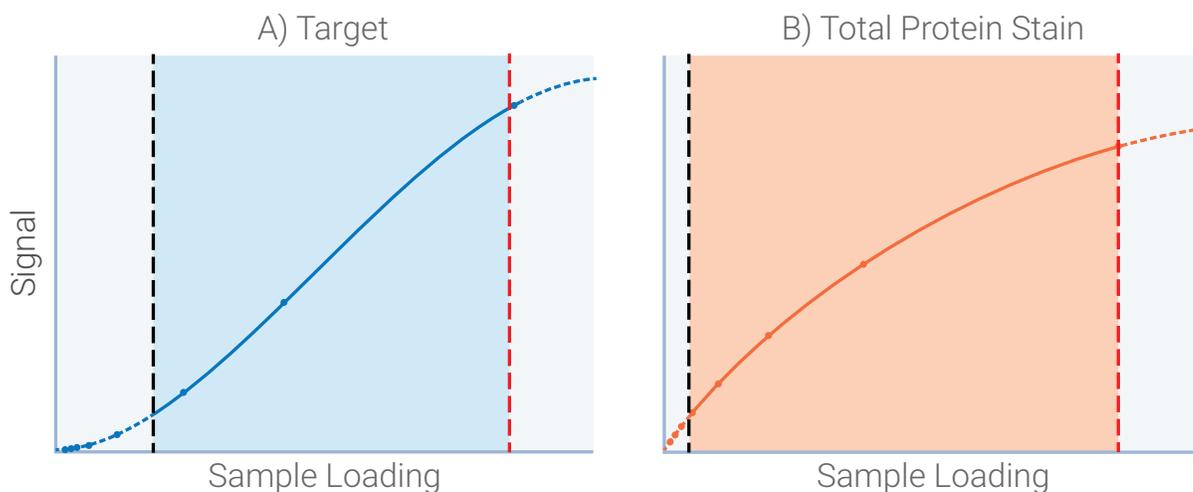


Figure 2. Estimate the boundaries for the linear range of sample loading. The estimated lower end (black dashed lines) and upper end (red dashed lines) are indicated for the target protein (A) and total protein stain (B). Shaded regions indicate the range of sample loading that produces a linear signal response on each graph. Linear range boundaries may be different for the target and total protein stain.

3. Combine the linear ranges of sample loading to identify the available range where the target and total protein stain can both be detected with a linear signal response (Figure 3). Use a secondary y-axis to plot the 700 nm and 800 nm signals on the same graph.
4. For your experiments, choose a level of sample loading that falls in the middle of the combined linear range for the target protein and Revert™ 700 Total Protein Stain (X, Figure 3). Avoid working near the boundaries of the combined range.

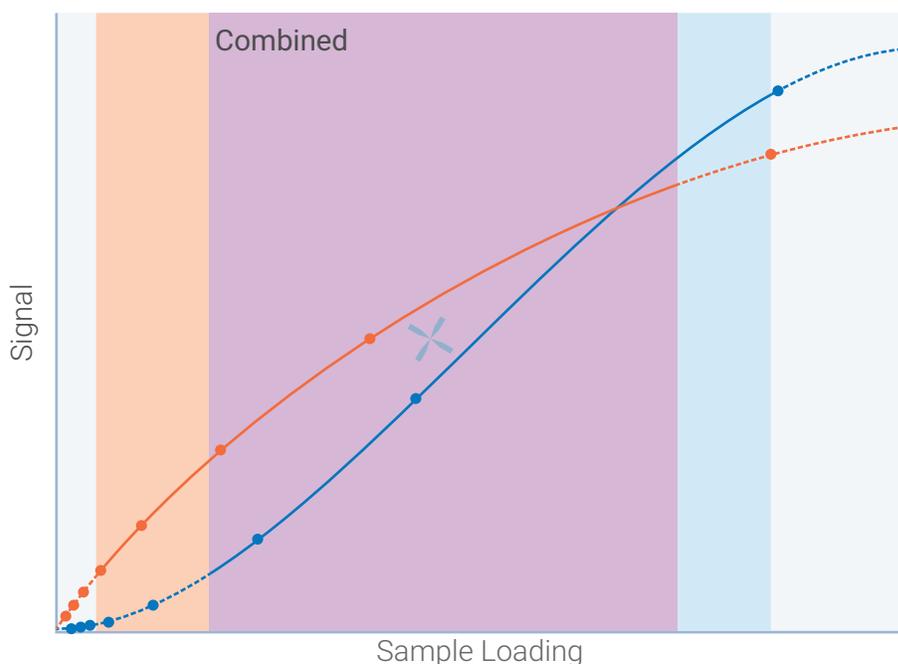


Figure 3. Determine the appropriate amount of sample to load. This combined graph indicates the appropriate range of sample loading for accurate detection of the target and total protein stain in a QWB experiment (X, center).

Note: If the target protein is strongly upregulated or downregulated, you should examine the linear range of sample loading for both untreated and treated samples.

VI. Protocol: Determining the Linear Range for a Target Protein and HKP

More Info: Empiria Studio® Software provides a dedicated workflow for this process. Go to licor.com/empiria to learn more.

Follow the instructions in this section if a housekeeping protein will be used as the internal loading control for QWB normalization. This section also applies to normalization with a pan-specific antibody for analysis of phosphorylation or other post-translational modifications. If you plan to use a total protein stain for normalization, see Section V.

1. Perform a Western blot with a serial dilution of sample (an 8-12 point, two-fold serial dilution is a good place to start).

Note: Make sure the dilution series includes sample loading amounts above and below the amount of sample you expect to load in each lane.

2. Prepare samples to be loaded on the gel with sample loading buffer.
3. Denature sample by heating at 95 °C for 3 min (or 70 °C for 10 min).
4. Load a uniform amount of sample protein in each lane.
5. Separate protein by SDS-PAGE.
6. Transfer proteins to immobilizing membrane.
7. Perform Western blot detection of target protein and HKP, according to the *Near-Infrared Western Blot Detection Protocol* (licor.com/NIRWesternProtocol; LI-COR).

Note: Use the 800 nm channel to detect the target protein. This channel provides the lowest membrane background and will maximize the sensitivity of detection.

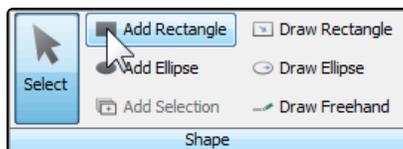
8. Image membrane with an Odyssey imaging system (Classic, CLx or Fc) in the 700 and 800 nm channels.
Adjust settings so that no saturation appears in the bands to be quantified.

Target Protein and HKP Quantification

Use Image Studio™ Software (licor.com/islite) to quantify the fluorescent signals of the HKP (700 nm) and target protein (800 nm).

HKP Quantification (700 nm)

1. Select and view the 700 nm channel image only.
2. Add shapes to bands.
 - a. In the Shape group on the Analysis tab, click Add Rectangle.



- b. Click each band to be analyzed, and an appropriately sized shape will be added around the band.

More Info: For help choosing the right background subtraction method, see licor.com/BgSubtractHelp.

3. Export the quantification data for your HKP.
 - a. Click Shapes to open the Shapes data table.

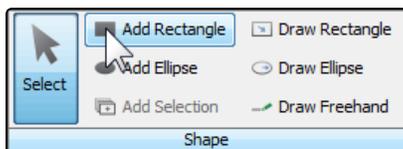


- b. Select shape data, then copy and paste data into a spreadsheet.

Note: All data fields will be exported, but "Signal" is the field of interest for analysis.

Target Protein Quantification (800 nm)

1. Select and view the 800 nm channel image only.
2. Add shapes to bands.
 - a. In the Shape group on the Analysis tab, click Add Rectangle.



- b. Click each band to be analyzed, and an appropriately sized shape will be added around the band.

More Info: For help choosing the right background subtraction method, see licor.com/BgSubtractHelp.

3. Export the quantification data for your target.
 - a. Click Shapes to open the Shapes data table.



- b. Select shape data, then copy and paste data into a spreadsheet.

Note: All data fields will be exported, but “Signal” is the field of interest for analysis.

Estimate Combined Linear Range for Target and HKP

1. Plot signal intensity vs. sample loading for the target and HKP (as shown in Figure 4).
2. Estimate the boundaries of the linear range of sample loading for the target and HKP (shaded regions, Figure 4).
 - a. Estimate the low end of the linear range by selecting a level of sample loading that is above the level of membrane background (black dashed lines, Figure 4).
 - b. Estimate the upper end of the linear range by selecting a level of sample loading that is below the region where band intensity begins to plateau due to saturation (red dashed lines, Figure 4).

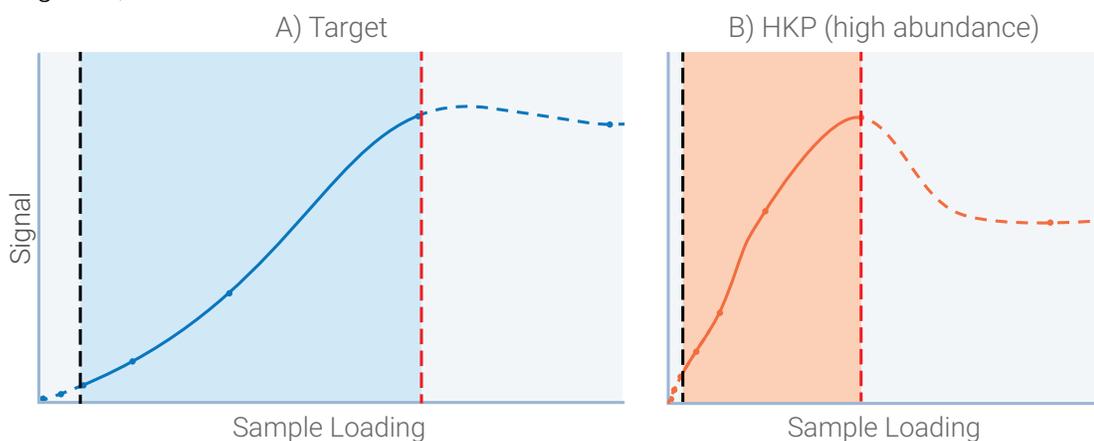


Figure 4. Estimate the boundaries for the linear range of sample loading. The estimated lower end (black dashed lines) and upper end (red dashed lines) are indicated for the target protein (A) and abundant HKP (B). Shaded regions indicate the range of sample loading that produces a linear signal response on each graph. Linear range boundaries may be very different for the target and HKP.

3. Combine the linear ranges of sample loading to identify the available range where the target and HKP can both be detected with a linear signal response (Figure 5). Use a secondary y-axis to plot the 700 nm and 800 nm signals on the same graph.
4. For your experiments, choose a level of sample loading that falls in the middle of the combined linear range for the target protein and HKP (X, Figure 5). Avoid working near the boundaries of the combined range.

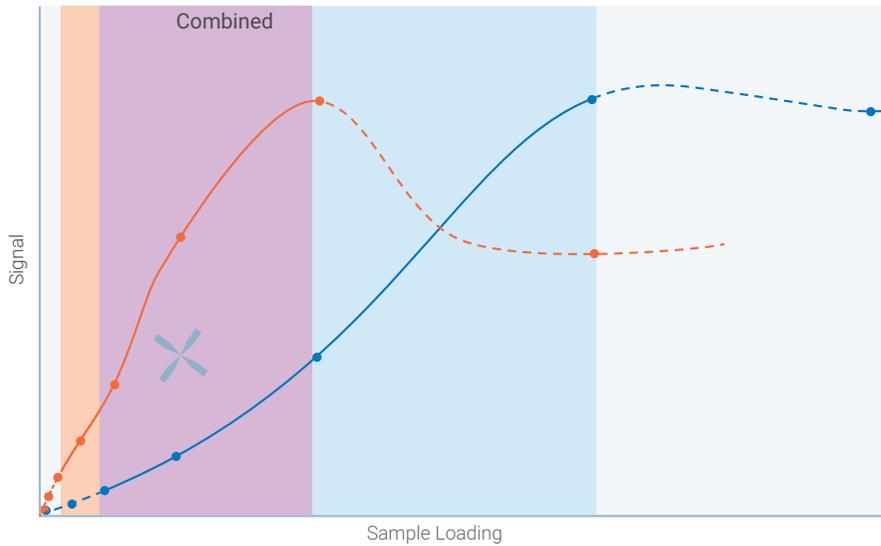


Figure 5. Determine the appropriate amount of sample to load. This combined graph indicates the appropriate range of sample loading for accurate detection of the target and HKP loading control in a QWB experiment (X, center).

5. If the target is low in abundance and the HKP is highly expressed, it may not be possible to identify a level of sample loading that is appropriate for both proteins. Consider using a total protein stain as an internal loading control for QWB normalization.

Note: If the target protein is strongly upregulated or downregulated, you should examine the linear range of sample loading for both untreated and treated samples.

VII. Further Reading

Please see the following for more information about QWB analysis.

Western Blot Normalization Handbook

licor.com/handbook

The *Normalization Handbook* describes how to choose and validate an appropriate internal loading control for normalization.

Good Normalization Gone Bad

licor.com/GNGB

Good Normalization Gone Bad presents examples of normalization that have been adversely affected by common pitfalls and offers potential solutions.

Western Blot Normalization White Paper

licor.com/normalizationreview

This white paper comprehensively reviews the literature of Western blot normalization.

Determining the Linear Range for Quantitative Western Blot Detection

licor.com/LinearRange

This protocol explains how to choose an appropriate amount of sample to load for QWB analysis.

Revert™ 700 Total Protein Stain Normalization Protocol

licor.com/RevertNormalization

This protocol describes how to use Revert 700 Total Protein Stain for Western blot normalization and analysis.

Pan/Phospho Analysis For Western Blot Normalization

licor.com/PanProteinNormalization

This protocol describes how to use pan-specific antibodies as an internal loading control for normalization.

Housekeeping Protein Validation Protocol

licor.com/HKP-Validation

This protocol explains how to validate an HKP for use as an internal loading control, by demonstrating that HKP expression is stable in the relevant experimental samples.

Housekeeping Protein Normalization Protocol

licor.com/HKP-Normalization

This protocol describes how to use a housekeeping protein for Western blot normalization and quantitative analysis.

References

1. Robasky, K, Lewis NE, and Church GM. *Nat. Rev. Genet.* 15: 56–62 (2014).
<http://www.nature.com/nmeth/journal/v11/n9/pdf/nmeth.3091.pdf>
2. Naegle K, Gough NR, and Yaffe MB. *Sci Signal.* 8:fs7 (2015).
<https://www.ncbi.nlm.nih.gov/pubmed/25852186>



© 2019 LI-COR, Inc. LI-COR, Odyssey, Revert, Empiria Studio, and Image Studio are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH

Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.

St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com