

# Establishing Image Integrity in Western Blotting: A Practical Guide to Documentation and Data Verification

Technical Overview

Nikolas Chmiel and John Riggs

Bio-Rad Laboratories, Inc.

Western blotting remains one of the most widely used techniques for protein analysis and detection, despite advances in mass spectrometry and other proteomics methods. However, image integrity and proper documentation of western blots have become increasingly critical for publication and scientific reproducibility. Major journals, funding agencies, and institutional review bodies now require rigorous evidence that blot images have not been manipulated or misrepresented. This article discusses practical approaches to establishing, documenting, and verifying image integrity in western blotting experiments, including free and accessible tools available to all researchers.

The scientific community has identified several high-profile cases of image manipulation in published western blots, ranging from simple copy-paste artifacts to more subtle alterations of contrast and brightness that misrepresent biological findings. These incidents have prompted journals including *Nature*, *Cell*, the *Journal of Biological Chemistry*, and the *Journal of Cell Biology* to implement mandatory image integrity checks during peer review. Establishing robust documentation practices is now essential for any researcher seeking to publish gel-based data.

## Understanding Image Integrity in Western Blotting: What Constitutes Image Manipulation?

Image integrity violations in western blots can fall into several categories:

- Cropping images to remove bands that indicate inconvenient results
- Selectively removing or duplicating lanes or lane portions
- Applying inconsistent brightness and contrast adjustments, even minor ones, leads to improper band assignment and misrepresents the relative intensity of bands, which is the quantitative foundation of western blot analysis
- Removing unwanted data with blending or cloning tools
- Combining lanes from different gels or exposures without clear documentation
- Other alterations of band appearance, such as rotating bands

Importantly, many image manipulation practices are performed inadvertently. Non-destructive adjustments such as exposure time optimization, background subtraction, and image rotation for presentation purposes are scientifically reasonable and necessary; however, these must be applied uniformly across all lanes being compared and must be fully disclosed in figure legends or methods.

## Documentation as Evidence of Data Integrity

The most practical and defensible approach to demonstrating image integrity is comprehensive documentation. This includes:

- The original, unprocessed image (usually the TIFF file generated directly by the imaging system)
- The processed image(s) submitted for publication with clear notation of any adjustments applied
- The specific software and parameters used for any adjustments
- Evidence that identical adjustments were applied to all samples being quantitatively compared

Journal editors and peer reviewers increasingly request these source files, either in supplementary materials or via the author's institutional repository. Having complete documentation readily available expedites the review process, strengthens the credibility of the work, and protects the researcher from questions about data integrity.



## Practical Documentation Workflow

### Step 1: Acquire and Archive the Original Image

Most gel imaging systems (Bio-Rad, LiCor, Cytiva) save images in proprietary formats that can be exported to lossless formats such as TIFF. Immediately after image acquisition, export the raw image as an uncompressed TIFF without any gain or contrast adjustments. Store this file in a laboratory archive with a clear naming convention (e.g., *20260515\_WB\_p53\_exp1\_lane1-6\_raw.tif*). This becomes the ultimate reference document. Do not open or edit this file after archiving; instead, work from a copy. The copy, in its native, proprietary format, can be used for further analysis in tools such as [Image Lab Software](#).

**Note:** *This workflow assumes access to institutional storage infrastructure and imaging systems under direct investigator control. Researchers using shared core facilities may have limited access to raw proprietary files or export settings; in such cases, work closely with core facility staff to establish archiving protocols and request maximum resolution exports in standardized formats.*

### Step 2: Document Acquisition Parameters

Create a simple text file or spreadsheet entry to record: exposure time, gain setting, laser power (if applicable), filter settings, and bit depth. For example, *ChemiDoc System imaging: 1 sec exposure, 1.0 gain, auto exposure off, 16-bit*. These metadata are essential for understanding the quantitative relationship between pixel intensity and protein amount. Bit depth is particularly critical. 16-bit images provide a greater dynamic range than 8-bit images and are preferred for quantitative analysis. Of note, most modern imaging systems are 16-bit imagers.

**Note:** *For some western blot imagers, some of the aforementioned settings may not be user-controlled and may be managed directly by the system. Depending on the imaging system used, the metadata information may be available and automatically stored within the proprietary image format.*

### Step 3: Apply Uniform Adjustments for Presentation

For figures intended for publication, any brightness or contrast adjustments should be documented and applied uniformly to all lanes being quantitatively compared. Many journals accept images processed with standard software (ImageJ, Python with PIL/scikit-image) provided adjustments are clearly reported. Document the specific adjustment. Example: Brightness +5, Contrast +10 applied uniformly to all lanes.

**Caveat:** *"Uniform" adjustments can present practical challenges in real blots where background fluorescence varies across the gel due to uneven transfer or edge effects. In such cases, researchers may find that a single adjustment setting does not optimally visualize all bands. Best practice is to perform all quantitative densitometry analysis on the raw, unadjusted image; adjustments are acceptable only for publication clarity.*

### Step 4: Create and Label the Publication Figure

The figure submitted for publication should include a clear legend:

- Antibody, including vendor, catalog, and clone number
- Band sizes and molecular weight markers
- Any image processing applied. Example: Uniform brightness and contrast adjustment applied to all lanes for clarity
- The normalization method used (housekeeping protein, total protein, or both). Example: western blot probed with anti-p53 (DO-1, 1:1000) and stripped/reprobed with anti-GAPDH (14C10, 1:3000) as loading control

## Quantification and Total Protein Normalization

Even with careful documentation and transparent image handling, the reliability of western blot data ultimately depends on how those images are quantified and interpreted. Normalization is therefore a critical extension of image integrity, ensuring that observed differences reflect biology rather than variability in sample loading or analysis.

In practice, this has traditionally been achieved through densitometry, with band intensity normalized to a single loading control, typically a housekeeping protein such as GAPDH or actin.

This approach remains widely used but has recognized limitations. Single-control normalization assumes that the loading control protein is unchanged across experimental conditions, and that total protein loading is consistent across lanes. These assumptions require validation for each experimental system.

### Total Protein Normalization and Total Protein Loading (Stain-Free and Related Methods)

An alternative normalization approach uses total protein detection via Stain-Free technology (e.g., Bio-Rad **Stain-Free** gels) or reversible staining methods such as Ponceau S. The Stain-Free technology approach exploits the intrinsic fluorescence of tryptophan residues in proteins under UV activation, allowing detection and quantification of total protein per lane without additional staining. The stated advantages include:

- Normalization to total protein rather than a single control
- Detection of potential loading inconsistencies that might be missed with a single housekeeping protein
- Reduction in membrane stripping and reprobing steps

Recent evidence supports improved reproducibility with total protein normalization. An analysis comparing normalization methods (Maloy et al, 2023) found that Stain-Free normalization reduced variability compared to actin or  $\beta$ -tubulin normalization and reduced the sample size needed for statistical significance by >50%.

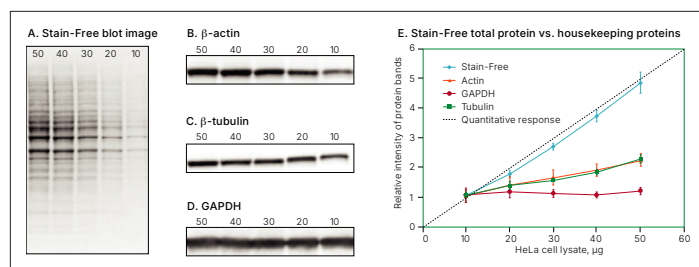
**Note:** *Stain-Free methodology offers a broad dynamic range and typically aligns well with typical protein load and reduced saturation issues commonly found with housekeeping proteins. Stain-Free total protein normalization, like any other normalization technique, needs to be validated by the user.*

### Important Caveats Regarding Total Protein Normalization

- Stain-Free fluorescence efficiency varies by protein context. Aromatic residue accessibility, hydrophobic environment, and neighboring charges all affect tryptophan fluorescence; equal concentrations of different proteins do not necessarily produce equal fluorescent signals
- Stain-Free fluorescence is optimized for samples containing ~10  $\mu$ g or more of protein
- Pretransfer Stain-Free images measure total protein on the gel before transfer; post-transfer methods assume equal transfer efficiency across lanes, which is not guaranteed if some lanes contain unusual protein mixtures or high salt. The Stain-Free workflow addresses this by enabling real-time quality control of the western blotting process and capturing a post-transfer total protein image of the blot for normalization
- Global protein changes can erase biological significance. If your experimental treatment alters total protein abundance (e.g., protein synthesis inhibition, starvation, differentiation), normalizing to total protein can obscure or reduce the apparent effect. In such cases, housekeeping protein normalization or dual normalization (both total protein and housekeeping control) may be more appropriate
- Standardization across platforms remains incomplete. Different manufacturers, gel types, membrane types, and imaging systems give different Stain-Free signal relationships to protein amount. The method is not yet standardized across all equipment combinations

For publication, the *Journal of Biological Chemistry* strongly recommends normalization to total protein but emphasizes that housekeeping proteins should be used only if the investigator has demonstrated that their expression is unaffected by experimental treatments. The most rigorous approach is reporting both total protein normalized and housekeeping protein normalized values in supplementary data, allowing readers to assess the robustness of conclusions.

For more information on journal requirements for publication, visit <https://www.bioradiations.com/from-gel-to-journal-a-practical-guide-to-western-blot-submission/Radiations>



**Fig. 1. Stain-Free technology provides a 1:1 correspondence between total protein and signal over a wide range of protein concentrations.** **A**, Stain-Free total protein signal; **B**,  $\beta$ -actin signal; **C**,  $\beta$ -tubulin signal; **D**, GAPDH signal; **E**, comparison of linearity of total protein normalization vs. housekeeping proteins over a range of protein concentrations.

## Freely Available Tools for Image Analysis and Documentation

### ImageJ (NIH)

ImageJ ([imagej.net](http://imagej.net)) is a public-domain image analysis software developed by the National Institutes of Health. It is widely accepted by major journals for gel image analysis and quantification. Key capabilities relevant to western blotting include:

Feature	Description
Densitometry	Quantify band intensity and area; generate pixel intensity profiles
ROI Selection	Define regions of interest (bands) with precision; automated lane detection available
Background Subtraction	Remove background fluorescence with rolling-ball or other algorithms
Documentation & Export	Save analysis results, overlay selections on images, export high-resolution figures

ImageJ is particularly valuable because it preserves full metadata and allows creation of macros to apply identical processing to multiple images. However, users should be aware that ImageJ has a learning curve, and macro consistency depends on proper validation against known standards. Fiji, an ImageJ distribution ([fiji.sc](http://fiji.sc)), includes additional plugins and is recommended for new users seeking a more integrated environment.

### Other Freely Available Resources

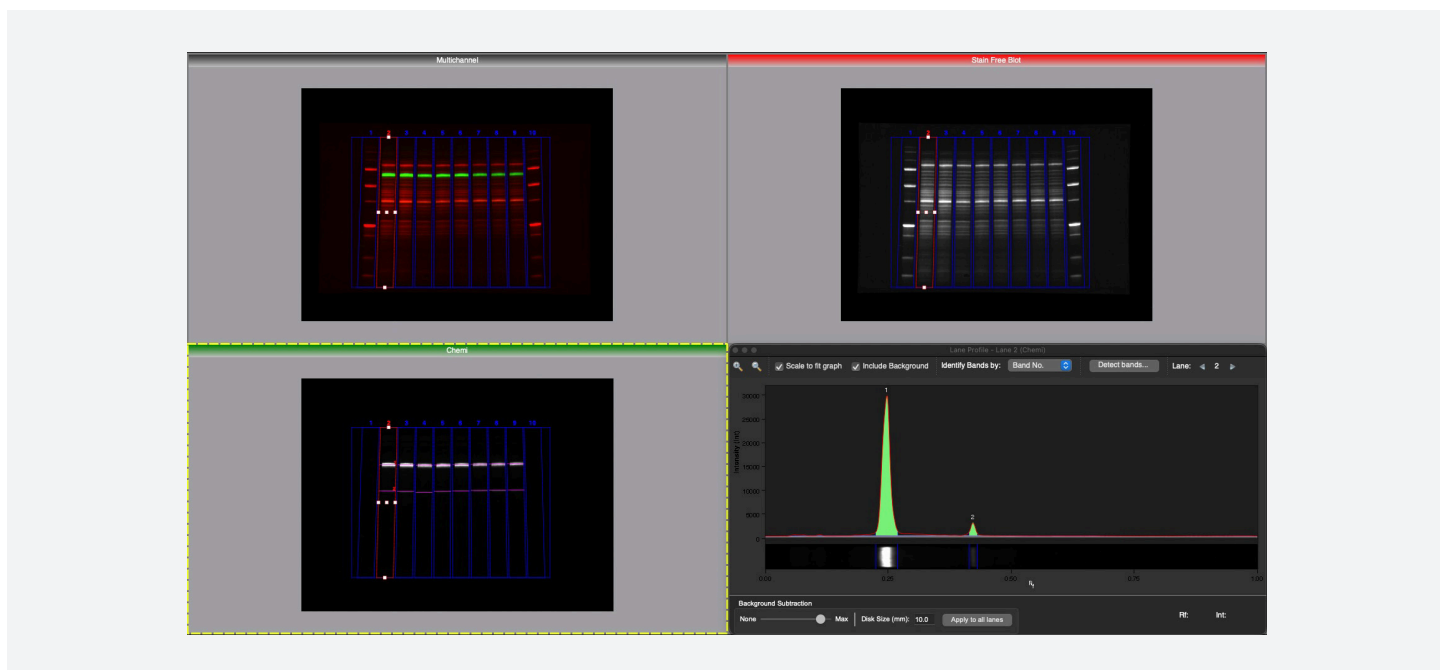
**GIMP (GNU Image Manipulation Program):** general-purpose image editor ([gimp.org](http://gimp.org)) useful for cropping, documenting regions, and adding annotations to gel images. Not specifically designed for scientific quantification, but appropriate for documentation and figure preparation.

**Python with scikit-image:** for researchers comfortable with programming, Python libraries (scikit-image, OpenCV, scipy) provide full control over image processing pipelines. Scripts can be version-controlled and shared, enhancing reproducibility. However, Python-based analysis requires coding expertise and adequate validation.

**R and EBImage:** the Bioconductor project provides EBImage and related packages for gel image analysis within the R statistical environment, allowing seamless integration with downstream statistical analysis.

### Bio-Rad Software Solutions

**Image Lab Software** supports image integrity by preserving the full context of data from acquisition through analysis within a secure, traceable environment. By maintaining native file formats that embed acquisition metadata and analysis parameters, the software enables users to document exactly how images were generated and processed, ensuring that any adjustments are transparent and reproducible. Features such as controlled export to lossless formats, audit-ready records, and consistent application of analysis settings across samples help researchers meet increasing expectations for documentation, reinforcing confidence that reported results faithfully reflect the original data.



## Practical Checklist for Submitting Western Blot Data

Before submitting a manuscript containing western blot data, use the following checklist to ensure image documentation and reporting are complete and transparent:

- Archive original (raw) image files from the imager and the lossless format (TIFF) with clear naming and acquisition metadata
- Record the acquisition parameters: exposure time, gain, laser power, filter settings, bit depth, membrane type, and binning
- Document the normalization method and rationale: total protein, housekeeping protein, or both
- Clearly note any image processing applied (software, specific parameters) and ensure it has been applied consistently across compared lanes
- Ensure that figure legends include antibody names and sources, including vendor name, clones, and catalog numbers. Also include band sizes, molecular weight markers, and normalization method
- Thoroughly describe the western blot protocol, antibody dilutions, detection method, quantification approach, and number of biological/technical replicates in the methods section
- Report densitometric analysis results (raw and normalized values) in supplementary tables, including all individual replicates
- Ensure that no lanes have been deleted or spliced between different blots without clear labeling and explicit description
- Include full blot images (uncropped) in supplementary materials
- If total protein normalization is used, include images and densitometric analysis of both total protein detection and target protein antibody probing
- Be ready to provide source files, analysis code/macros, and additional details upon reviewer request

### Limitations of the Checklist Approach

While comprehensive documentation is essential, it is important to recognize that passing a documentation checklist does not guarantee either biological validity or quantitative accuracy. The following limitations should be considered:

- Documentation is not QA/QC. Complete metadata, archived raw images, and transparent reporting satisfy publication standards but do not address whether the underlying experimental design, biological replicates, or statistical analysis are sound
- Normalization method choice is not neutral. Documentation of which normalization method you used does not prove you chose correctly for your system. Total protein normalization can obscure treatment-induced global protein changes; housekeeping protein normalization can miss those same changes. The checklist requires reporting the method, not validating it
- "Uniformly applied" adjustments do not address intrinsic gel heterogeneity. Uneven transfer, regional background differences, and edge effects can make truly uniform adjustments impractical. Best practice is performing densitometry on raw images, but the checklist may encourage researchers to optimize display at the expense of analytical rigor
- A study that passes the image integrity checklist is one piece of achieving overall scientific integrity. Scientists should still report additional experimental information including sample source, number of technical and biological replicates, preparation conditions, and other details
- Institutional constraints can limit archiving feasibility. Researchers using shared core facilities, limited storage systems, or equipment purchased over 10 years ago may lack access to lossless raw images or detailed acquisition parameters. The checklist assumes infrastructure many labs do not have
- Supplementary data are often not peer-reviewed. Supplementary raw images may be uploaded without verification that they match the main figures or that densitometry was performed correctly. Journals often do not screen supplementary files as rigorously as main text

In summary, the documentation checklist is a necessary but not sufficient condition for publication-quality western blot data. Researchers should view the checklist as a foundation for transparency, not a substitute for thoughtful experimental design, appropriate controls, and careful interpretation of results.

## Conclusion

Rigorous documentation of image integrity in western blotting is now expected by journals, funders, and the scientific community. The practical approach is straightforward: archive the original image, document all processing steps, apply adjustments consistently, and transparently report the methods and results. Researchers should understand, however, that documentation and transparency are necessary but not sufficient for scientific validity. Careful selection of normalization methods, consideration of biological context, and appropriate statistical analysis remain essential.

Recent evidence increasingly supports total protein normalization (via Stain-Free or reversible staining) over single housekeeping proteins as a more robust approach, with potential for reduced sample sizes and improved reproducibility. However, this advantage is context-dependent: total protein normalization may not be universally appropriate when the treatment itself changes global protein levels, and accuracy may vary depending on the total protein normalization methods applied. Researchers should validate their chosen normalization strategy for their specific biological system rather than assuming that any approach is universally correct.

Excellent free tools such as ImageJ are available for image analysis and quantification, eliminating cost as a barrier to rigorous work. Of note, Stain-Free total protein normalization using proprietary Stain-Free technology from Bio-Rad requires the use of Image Lab Software, which is included with every GelDoc™ Go Gel Imaging System\* and every ChemiDoc™ Imaging System.\* The time invested in establishing proper documentation and validation early in research, rather than scrambling to reconstruct methods after data collection, is the most efficient approach and strengthens both the quality and defensibility of published results.

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

Visit [bio-rad.com/GoStainFree](https://www.bio-rad.com/GoStainFree) for more information.

Bio-Rad, ChemiDoc, and GelDoc are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.  
© 2026 Bio-Rad Laboratories, Inc.

## Further Reading

- Cromey DW (2013). Digital images are data: and should be treated as such. *Methods Mol Biol* 931, 1-27.
- Eaton SL Roche et al. (2013). Total protein analysis as a reliable loading control for quantitative fluorescent western blotting. *PLoS One* 8, e72457.
- Fosang AJ and Colbran RJ. (2015). Transparency is the key to quality. *J Biol Chem* 290, 29692-29694.
- Gassmann M et al. (2009). Quantifying Western blots: pitfalls of densitometry. *Electrophoresis* 30, 1845-1855.
- Gilda JE and Gomes AV (2013). Stain-free total protein staining is a superior loading control to  $\beta$ -actin for Western blots. *Anal Biochem* 440, 186-188.
- Maloy A et al. (2023). Stain-Free total-protein normalization enhances the reproducibility of Western blot data." *Anal Biochem* 654, 114840.
- Rossner M and Yamada KM (2004). What's in a picture? The temptation of image manipulation. *J Cell Biol* 166, 11-15.
- Taylor SC (2013). A defined methodology for reliable quantification of Western blot data. *Mol Biotechnol* 55, 217-226.
- Wang Q et al. (2023). Western blot normalization: Time to choose a proper loading control seriously. *Electrophoresis* 44, (854-863).
- Westerberg LJS et al. (2025). Superior normalization using total protein for western blot analysis of human adipocytes. *PLoS One* 18, e0328136.

## Free Resources and Tools

- ImageJ/Fiji: [imagej.net](https://imagej.net) and [fiji.sc](https://fiji.sc) (Essential open-source image analysis software; Fiji recommended for new users.)
- GIMP: [gimp.org](https://gimp.org) (General-purpose image editor useful for documentation and annotation.)
- Bioconductor: [bioconductor.org](https://bioconductor.org) (Open-source R packages for computational biology, including image analysis integration.)
- scikit-image: [scikit-image.org](https://scikit-image.org) (Python library for image processing and analysis.)
- EMBO Resource Center Guidelines: [embopress.org](https://embopress.org) (EMBO Journal's detailed image integrity guidelines.)
- Journal of Biological Chemistry Image Guidelines: [jbc.org](https://jbc.org) (Comprehensive guidance on normalization methods and supporting data requirements.)



**Bio-Rad**  
Laboratories, Inc.

Life Science  
Group

Website [bio-rad.com](https://www.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 55 11 3065 7550 Canada 1 800 361 1808 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 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 000 800 00 24 67 23 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 55 5488 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 7 495 721 14 04 Singapore 65 6415-3170 South Africa 27 21 531 7504 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 662 651 8311 United Arab Emirates 971 4 818 7300 United Kingdom 00 800 00 24 67 23