

ReadiPrep™ Western Blot Stripping Buffer

Catalog number: 60513
Unit size: 100 mL

Component	Storage	Amount (Cat No. 60513)
ReadiPrep™ Western Blot Stripping Buffer	Refrigerated (2-8 °C), Minimize light exposure	100 mL

OVERVIEW

ReadiPrep™ Western Blot Stripping Buffer is a ready-to-use solution for removing primary and secondary antibodies from Western blot membranes without damaging immobilized proteins. This buffer enables efficient stripping of chemiluminescent, fluorescent, and chromogenic antibody complexes, allowing the same blot to be reprobed multiple times. Its mild formulation preserves protein integrity and membrane binding, making it ideal for applications requiring sequential detection of different targets. The protocol requires only a short room temperature incubation and is compatible with PVDF and nitrocellulose membranes. ReadiPrep™ Western Blot Stripping Buffer reduces reagent use and increases data output from a single blot.

AT A GLANCE

1. Wash blot with PBS or TBS for 5 minutes to remove substrates.
2. Incubate blot in ReadiPrep™ Western Blot Stripping Buffer for 20–60 minutes at room temperature (optional 37°C for tough antibodies).
3. Wash twice with PBS or TBS, then proceed with reprobing.

SAMPLE EXPERIMENTAL PROTOCOL

1. Wash the western blot for 5 minutes with PBS or TBS to remove the chemiluminescent substrates.
2. Add sufficient ReadiPrep™ Western Blot Stripping Buffer to cover the blot and incubate for 20 to 60 minutes at room temperature with gentle shaking.
Note: Optimization of both incubation time and temperature is essential for best results. For some antibodies, room temperature incubation is sufficient. However, high-affinity antibodies or saturated blots may require incubation for an additional 5 to 10 minutes at 37°C.
3. Remove the blot from the ReadiPrep™ Western Blot Stripping Buffer.
4. Wash blot 2 times with PBS or TBS for 5 minutes each with gentle shaking.
5. After the wash, blot can be reused for a second immunoprobing experiment.

Note: To determine the successful stripping of antibodies, HRP or AP labeled secondary antibody can be added followed by addition of new chemiluminescence substrates. If no signal is detected then the primary antibody has been successfully removed from the antigen.

IMPORTANT NOTES

1. ReadiPrep™ Western Blot Stripping Buffer is not recommended to be used for stripping fluorescent western blots. The buffer is compatible with most chemiluminescence substrates.
2. Reblocking the membrane is recommended after stripping
3. ReadiPrep™ Western Blot Stripping Buffer is supplied in ready-to-use format. There is no need to dilute the buffer.
4. For performing multiple strippings and reprobing, it is

- recommended to probe for low abundant proteins first.
5. Bring the ReadiPrep™ Western Blot Stripping Buffer to room temperature prior to use.

EXAMPLE DATA ANALYSIS AND FIGURES

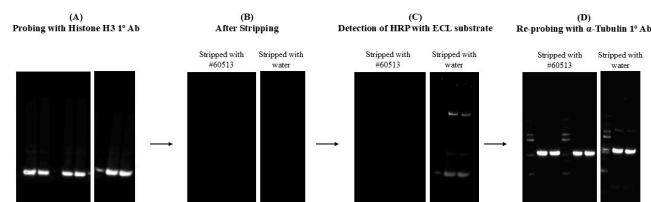


Figure 1. Untreated HeLa cell lysate (10 µg) was separated by SDS-PAGE and transferred to nitrocellulose membranes. Blots were probed with an Anti-Histone H3 antibody (Panel A). After stripping with ReadiPrep™ Western Blot Stripping Buffer (#60513), no residual signal was detected (Panel B). Signal was measured again after additional incubation of ECL substrates (Panel C). The blots were then re-probed with an Anti-Alpha-Tubulin antibody (Panel D). Amplitude® West ECL HRP Substrate (#26100) was used to develop the chemiluminescence response.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.