

Components

Components for the Quick Western Kit (926-69100):

- IRDye® 680RD Detection Reagent
- Intercept® (PBS) Blocking Buffer

Other Required Reagents

- 1X PBS
- Primary Antibody
- 1X PBS-T (1X PBS containing 0.1% Tween® 20)
- 20% Tween 20
- 20% SDS (If using PVDF membrane)

Specifications

- Fluorophore: IRDye 680RD
- Excitation Wavelength: 676 nm (in PBS)
- Emission Wavelength: 694 nm (in PBS)
- Form of Detection Reagent: IRDye 680RD Detection Reagent, lyophilized in water. Contains 10 mg BSA (IgG and protease free) per mg of Detection Reagent as a stabilizer and 0.01% sodium azide as a preservative.
- Storage: Store at 4 °C and protect from light. Will remain stable for up to 3 months at 4 °C in reconstituted form.
- Kit will perform approximately 25 Western blots.

Warning: *Contains sodium azide.*

Applications

The Quick Western Kit – IRDye 680RD provides a universal antibody detection reagent that can be combined with the primary antibody incubation step, eliminating the need for a secondary antibody. The overall time to complete a Western blot is reduced, while providing the advantages of near-infrared detection. The kit can be used to detect primary antibodies from a variety of hosts and has been shown to recognize primary antibodies to recombinant tagged proteins (e.g. 6X His, Myc, FLAG, etc.).

The kit also serves as a detection solution for post-immunoprecipitation samples by Western blot because it does not bind to denatured mouse monoclonal or rabbit monoclonal antibodies. The key benefit is the ability to use the same antibody for immunoprecipitation and post-immunoprecipitation detection by Western blot.

Note: *The IRDye 680RD Detection Reagent recognizes denatured polyclonal antibodies and is not recommended for detection of samples that have been immunoprecipitated using polyclonal antibody.*

Purity and Specificity

IRDye 680RD Detection Reagent is known to have high affinity for IgG from human, mouse, rabbit, guinea pig, goat, sheep, pig, cow, cat, dog, and donkey. The Detection Reagent is known to have lower affinity for mouse IgG₁, mouse IgA, rat, horse, and hamster IgGs and will not detect primary antibodies from a chicken host. The Detection Reagent has been specifically tested and qualified for Western blot applications. If additional affinity is required, please use IRDye conjugated secondary antibodies for detection.

Reconstitution

Reconstitute contents of vial with 0.25 mL 1X PBS. Mix gently by inverting. Protect from light and allow to rehydrate for at least 60 minutes before use. Centrifuge product if solution is not completely transparent after standing at room temperature.

Western Detection Protocol - Quick Western Kit

Step 1. Wet Membrane

For Nitrocellulose Membranes

Wet in 1X PBS for 5 minutes or until fully hydrated.

For Immobilon®-FL PVDF Membranes

1. Wet for 30 seconds in 100% methanol.
2. Wet in 1X PBS for 5 minutes.

Step 2. Block the Membrane

Place the membrane in an incubation box and block the membrane with Intercept Blocking Buffer for 1 hour at room temperature with gentle shaking.

Be sure to use sufficient blocking buffer to cover the membrane (a minimum of 0.4 mL/cm² is suggested).

Step 3. Prepare Detection Solution

1. Add primary antibody (dilute as recommended by vendor) to Intercept Blocking Buffer.
2. Add the IRDye 680RD Detection Reagent to the primary antibody + Intercept Blocking Buffer solution. Use 1 µL of IRDye 680RD to every 1 mL of Intercept Blocking Buffer (1:1000 dilution).
3. Add detergent:
 - For nitrocellulose membranes use 0.2% Tween® 20 in the detection solution.
 - For PVDF membranes use 0.2% Tween 20 and 0.02% SDS in the detection solution.

Note: Optimization is required to determine the final concentration of detergents. Concentration can range from 0.1 - 0.2% Tween 20 and 0.02 - 0.1% SDS.

Step 4. Decant and Discard Blocker

Step 5. Add Detection Solution

Add the detection solution prepared in Step 3.

Step 6. Incubate Blot in Detection Solution

Protect from light during incubation.

- Incubate using your standard primary antibody incubation time at room temperature on a platform shaker.
- If the standard primary antibody incubation time is overnight, store the blot at 4 °C on a platform shaker.

Step 7. Wash Membrane

1. Carefully decant detection solution.
2. Rinse the membrane in 15 mL of 1X PBS-T (1X PBS + 0.1% Tween 20).
3. Decant.
4. Add 15 mL of 1X PBS-T and incubate for 5 minutes at room temperature with gentle shaking. Protect from light.
5. Decant wash solution.
6. Repeat 2 additional times.

Step 8. Rinse Membrane

Rinse membrane with 15 mL of 1X PBS.

Step 9. Scan Membrane

Note: *Protect the membrane from light prior to scanning.*

Scan the membrane (wet or dry) on an Odyssey® Imaging System using the 700 nm channel.

Example

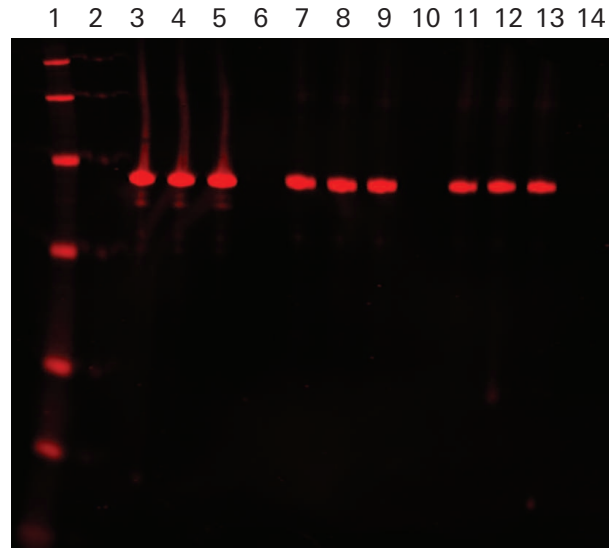


Figure 1. Primary antibody targeting β -actin was detected using IRDye[®] 680RD Detection reagent (1:1000; LI-COR P/N 926-69100) for 1 hour at room temperature. 5 μ g of A-10 (sc-3806), C32 (sc-2205), and HeLa (sc-2200) lysates were loaded on a 10% SDS-PAGE gel (NP0301BOX), in lanes 3-5, 7-9, and 11-13. Chameleon[®] Duo ladder (LI-COR P/N 928-70000) was loaded in lane 1. Lanes 2, 6, 10, and 14 are empty. The nitrocellulose membrane (LI-COR P/N 926-31090; 0.22 μ M; 7 cm x 8.5 cm) was blocked with Intercept (PBS) Blocking Buffer and β -actin (~45 kDa) was identified using Rabbit Anti- β -actin rabbit polyclonal antibody (LI-COR P/N 926-42210).

Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at licor.com/packinserts under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2021 LI-COR, Inc. LI-COR, Odyssey, Chameleon, Intercept, and IRDye 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