

## Protocol: Western Blot Protocol for p62/SQSMT1 Antibody (NBP1-48320)

### Materials

- 1X PBS
- Sample buffer, 2X Laemmli buffer: 4% SDS, 5% 2-mercaptoethanol (BME), 20% glycerol, 0.004% bromophenol blue, 0.125 M Tris HCl, pH 6.8
- 1X Running Buffer: 25 mM Tris-base, 192 mM glycine, 0.1% SDS. Adjust to pH 8.3
- 1X Transfer buffer (wet): 25 mM Tris-base, 192 mM glycine, 20% methanol Adjust to pH 8.3
- TBS
- TBST, TBS and 0.1% Tween
- Blocking solution: TBST, 5% non-fat dry milk
- rabbit anti-p62/SQSMT1 primary antibody (NBP1-48320) in blocking buffer (~2 µg/mL)

### Methods

1. Grow cells (e.g. HeLa or Neuro2A) *in vitro* to semi-confluency (70-75%).
2. Rinse cells with ice-cold 1X PBS and lyse cells with sample buffer.
3. Sonicate and incubate cells for 5 minutes at 95°C.

Tip: Cells are lysed directly in sample buffer.

Note: For cell lysis, an SDS containing buffer is recommended to identify the entire cellular pool of p62/SQSMT1.

4. Load 10-40 µg/lane of sample on a 12% polyacrylamide gel (SDS-PAGE).

Note: To determine autophagic flux based on p62/SQSTM1, immunoblot analysis should include samples treated with autophagy inducers and inhibitors.

5. Transfer proteins to a Nitrocellulose membrane for 60 minutes at 100V.

Tip: For more information on Western Blotting, see our Western Blot handbook: [https://images.novusbio.com/design/BR\\_westernblotguide\\_042816b.pdf](https://images.novusbio.com/design/BR_westernblotguide_042816b.pdf)

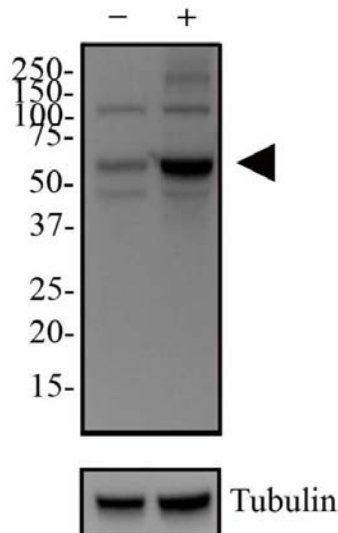
6. After transfer, rinse the membrane with dH<sub>2</sub>O and stain with Ponceau S for 1-2 minutes to confirm efficiency of protein transfer.
7. Rinse the membrane in dH<sub>2</sub>O to remove excess stain and mark the loaded lanes and molecular weight markers using a pencil.

8. Block the membrane using blocking buffer solution (5% BSA in TBST) for 1 hour at room temperature.
9. Rinse the membrane with TBST for 5 minutes.
10. Dilute the rabbit anti-p62/SQSTM1 primary antibody (NBP1-48320) in blocking buffer (~2  $\mu\text{g}/\text{mL}$ ) and incubate the membrane for 1 hour at room temperature.
11. Rinse the membrane with  $\text{dH}_2\text{O}$ .
12. Rinse the membrane with TBST, 3 times for 10 minutes each.
13. Incubate the membrane with diluted secondary antibody, according with product's specification, (e.g. anti-rabbit-IgG HRP-conjugated) in blocking buffer for 1 hour at room temperature.

Note: Tween-20 may be added to the blocking or antibody dilution buffer at a final concentration of 0.05-0.2%, provided it does not interfere with antibody-antigen binding.

14. Rinse the membrane with TBST, 3 times for 10 minutes each.
15. Apply the detection reagent of choice (e.g. BioFX Super Plus ECL) in accordance with the manufacturer's instructions.
16. Image the blot.

Note: p62/SQSTM1 is primarily degraded by autophagy, therefore its presence and amount in autophagosomes inversely correlates with autophagic activity. Inhibition of autophagosome content degradation increases the levels of p62/SQSTM1 detected by immunoblot.



*Western Blot: p62/SQSTM1 Antibody [NBP1-48320] - Cultured HeLa cells were treated with or without 50  $\mu$ M chloriquine for 24 hours as indicated. Cell lysates were prepared and separated on a 12% gel by SDS-PAGE. Protein was transferred to PVDF membrane and blocked in 5% non-fat milk. The membrane was then probed with 1  $\mu$ g/mL anti-p62/SQSMT1 in 1% milk and detected with an anti-rabbit HRP secondary antibody using chemiluminescence. Note the upregulation of p62 (arrowhead) in response to chloroquine treatment and the blockage of autophagy.*