

NB100-307 Protocol

Western Blot Protocol for ATM Antibody (NB100-307)

- 1) Resolve protein samples on a 6% SDS-PAGE gel at 185V for ~1.5 hours.
- 2) Transfer to PVDF membranes at 25V for ~1.5 hours.
- 3) Block the membrane with TBST+BSA and goat serum for 1 hour at RT.
- 4) Dilute purified primary antibody (NB100-307) to 1:1000 in blocking buffer.
- 5) Incubate membrane overnight at 4 degrees Celcius in diluted purified anti-ATM.
- 6) Wash 3 times ten minutes on a shaker.
- 7) Incubate membranes with HRP conjugated anti-mouse IgG for 1 hour (RT), diluted in blocking buffer.
- 8) Wash 3 times ten minutes on a shaker.
- 9) Add ECL reagent, as per kit directions, and expose for 1-5 seconds.

*The above information is only intended as a guide. The researcher should determine what protocol best meets their needs. Please follow safe laboratory procedures.