

NB100-306 Protocol

Western Blot Protocol for ATM Antibody (NB100-306)

Western Blot Procedure

- 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 primary antibody (NB 100-306) to 1:1,000 in blocking buffer.
- 5) Incubate membrane overnight at 4 degrees Celcius in diluted anti-ATM-kinase.
- 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.