

## NBP2-42215 Protocol

# Western Blot Protocol for BDNF Antibody (NBP2-42215)

### Reagents needed:

- Washing Buffer: Tris Buffer Saline with 0.01% of tween 20).
- Blocking Buffer: 5% skimmed milk powder in washing buffer).
- Secondary antibody, Horseradish peroxidase conjugated.
- Chemiluminescent solution (SuperSignal WestPico™, Pierce).

### Western blot Method:

- Perform SDS-PAGE using PVDF membrane. Cut into strips.
- Activate strips with methanol by dipping them into methanol for 5 min.
- Discard the methanol and take fresh methanol to repeat step b.
- Let the strips dry, and then add blocking solution and incubate at RT in a shaker for 30-45 minutes.
- Dilute primary antibody in blocking buffer. Incubate the number of strips required with the diluted primary antibody at room temperature for 2 hours in a shaker.
- Wash strips two times with washing buffer at 30 minutes intervals.
- Dilute HRP conjugated secondary antibody in blocking buffer. Add diluted secondary antibody to the membrane strips and incubate for exactly 1 hour while shaking at RT.
- Wash the strips with washing buffer for 2-3 hours with 3 to 4 changes on a shaker. This helps in reducing the back ground staining.
- Prepare the chemiluminescent solution (SuperSignal WestPico™) by mixing solution A and Solution B at 1:1. Mix well. Soak the strip in the chemiluminescent solution; keep for 3-5 minutes under constant shaking.
- Expose the membrane to a sheet of film and develop.