

NB100-105 Protocol

Immunohistochemistry Procedure (NB100-105)

Immunohistochemistry Procedure

1. If not previously done, bake sections at 60C for 30 minutes.
2. Hydrate sections through the following series.
 - A. 3 X 5 minutes xylenes
 - B. 3 X 5 minutes 100% EtOH
 - C. 2 minutes 95% EtOH
 - D. 2 minutes 70% EtOH
 - E. 1 minute 50% EtOH
 - F. 1 minute ddH₂O
 - G. 1 minute TBS
3. If dry sections are needed, circle sample with wax pencil.
4. Antigen unmasking was performed by microwaving in 0.1M sodium citrate (pH 6.0) for 2 X 5 minutes at power level 7. Cool for 15 minutes.
5. Rinse slides with TBS.
6. Quench slides in 0.3% hydrogen peroxide in MeOH (0.5 ml 30% stock in 50 ml MeOH) for 25 minutes.
7. Wash 2 X 5 minutes with TBS.
8. Block sections with 10% serum (from the host species of the secondary antibody) in TBS for 30 minutes.
9. Incubate NB 100-105 with sections overnight at 4C at a 1:50 dilution in 10% goat serum.
10. The following day, allow sections to sit at RT for 30 minutes.
11. Wash sections 3 X 5 minutes in TBS, followed by incubation with the secondary antibody diluted 1:100 in 10% goat serum for 30 minutes at RT.
12. After 15 minutes of incubation, make up ABC solution and allow to sit for 30 minutes.
13. Wash slides 3 X 5 minutes in TBS.
14. Block with ABC solution in TBS.
15. Make DAB solution.
16. Incubate with fresh DAB solution until signal develops, then place in ddH₂O. Dehydrate sections through graded alcohols to xylenes (reverse hydration steps) and coverslip with permount.