

NBP1-79010 Protocol

Immunohistochemistry-Paraffin protocol for CD117/c-kit Antibody (NBP1-79010)

IHC-P Protocol (NBP1-79010): https://www.novusbio.com/products/cd117-c-kit-antibody-r21-v_nbp1-79010

1. Deparaffinize the section in 3 changes of xylene, 5 minutes each.
2. Wash the section in 96%, 80% and 70% benzyl alcohol for 5 minutes each.
3. Rinse in distilled water.
4. Block the endogenous peroxidase by incubating the tissue in 3% hydrogen peroxide (H₂O₂) for 10 minutes.
5. Rinse in distilled water.
6. For antigen retrieval: immerse the slide in the citrate buffer, pH 6.0, 0.05% Tween-20*, and incubate in microwave (600W) for 10 minutes. (Alternatively adjust to your own protocol, keeping the required pH)
7. Remove the staining to room temperature and let the slide to cool (in citrate buffer, pH 6.0) for 15 minutes.
8. Rinse in distilled water.
9. Wash in 0.05 M Tris-HCl , pH 7.6 buffer supplemented with 0.2% of Tween-20 (buffer A) for 5 minutes.
10. Incubate the section with primary antibody in buffer A at the dilution 1:100 - 200 for 1 hour in the closed wet chamber.
11. Wash twice 5 minutes with buffer A.
12. Apply the secondary antibody (the protocol depends on the supplier), and proceed to standard immunohistochemistry protocol (HRP - Peroxide - DAB).
13. Wash twice 5 minutes with buffer A.
14. Apply the chromogen (DAB), 10 minutes.
15. Rinse in water.
16. Stain in hematoxylin for 5 minutes.
17. Wash in water - 10 minutes.
18. Dehydrate the section in 2 changes of 96% benzyl alcohol for 5 minutes each.
19. Wash the section in 2 changes of xylene for 2 minutes each.
20. Mount the slide for observation.