

NLS5377 Protocol

Immunohistochemistry Protocol for Steroidogenic Factor 1 Antibody (NLS5377)

Immunohistochemistry Protocol for Steroidogenic Factor 1 Antibody (NLS5377):

https://www.novusbio.com/products/sf-1-nr5a1-steroidogenic-factor-1-antibody_nls5377

Immunohistochemistry

1. Prepare tissue with formalin fixation and by embedding it in paraffin wax.
 2. Make 4-mm sections and place on pre-cleaned and charged microscope slides.
 3. Heat in a tissue-drying oven - 45 minutes @ 60 degrees Celcius.
 4. Deparaffinize the tissues by wash drying the slides in 3 changes of xylene?? - 5 minutes each @ RT.
 5. Rehydrate the tissues by washing the slides in 3 changes of 100% alcohol for - 3 minutes each @ RT.
 6. Wash the slides in 2 changes of 95% alcohol for - 3 minutes each @ RT.
 7. Wash the slides in 1 change of 80% alcohol for - 3 minutes @ RT.
 8. Rinse the slides in gentle running distilled water - 5 minutes @ RT.
 9. Perform antigen retrieval by steaming the slides in 0.01M sodium citrate buffer (pH 6.0) - 20 minutes @ 99-100 degrees Celcius.
 10. Remove the slides from the heat and let stand in buffer - 20 minutes @ RT.
 11. Rinse the slides in 1X TBS-T - 1 minute @ RT.
- **Do not allow the tissues to dry at any time during the staining procedure****
12. Begin the immunostaining by applying a universal protein block -?? 20 minutes @ RT.
 13. Drain protein block from the slides and apply the diluted primary antibody -?? 45 minutes @ RT.
 14. Rinse the slide in 1X TBS-T -?? 1 minute @ RT.
 15. Apply a biotinylated anti-rabbit IgG (H+L) secondary?? - 30 minutes @ RT.
 16. Rinse the slide in 1X TBS-T -?? 1 minute @ RT.
 17. Apply an alkaline phosphatase steptavidin -?? 30 minutes @ RT.
 18. Rinse the slide in 1X TBS-T -?? 1 minute @ RT.
 19. Apply an alkaline phosphatase chromagen substrate -?? 30 minutes @ RT.
 20. Rinse the slide in distilled water - 1 minute @ RT.
- **This method should only be used if the chromagen substrate is alcohol insoluble (ie: Vector Red, DAB)****
21. Dehydrate the tissue by washing the slides in 2 changes of 80% alcohol for 1 minute each @ RT.
 22. Wash the slides in 2 changes of 95% alcohol for 1 minute each @ RT.
 23. Wash the slides in 3 changes of 100% alcohol for - 1 minute each @ RT.
 24. Wash the slides in 3 changes of xylene -?? 1 minute each @ RT.
 25. Apply cover slip.

"