

## NB100-2527 Protocol

### Immunohistochemistry-paraffin embedded sections

LOX Antibody: [https://www.novusbio.com/products/lox-antibody\\_nb100-2527](https://www.novusbio.com/products/lox-antibody_nb100-2527)

Antigen unmasking

1. Bring slides to a boil in 10 mM sodium citrate buffer pH 6.0 then maintain at a sub-boiling temperature for 10 minutes.
2. Cool slides on bench top for 30 minutes.

#### Staining

1. Wash sections in dH<sub>2</sub>O three times for 5 minutes each.
2. Wash section in wash buffer (1X PBS/0.1% Tween-20 (1X PBST)) for 5 minutes.
3. Block each section with 100-400ul blocking solution (1X PBST, 5% goat serum) for 1 hour at room temperature.
4. Remove blocking solution and add 100-400 ul primary antibody diluted in 1X PBST, 5% goat serum to each section.
5. Incubate overnight at 4C.
6. Remove antibody solution and wash sections in wash buffer three times for 5 minutes each.
7. Add 100-400 ul biotinylated secondary antibody, diluted in 1X PBST, 5% goat serum.
8. Incubate 30 minutes at room temperature.
9. Remove secondary antibody solution and wash sections three times with wash buffer for 5 minutes each.
10. Add 100-400 ul Streptavidin HRP reagent to each section and incubate for 30 minutes at room temperature.
11. Wash sections three times in wash buffer for 5 minutes each.
12. Add 100-400 ul DAB substrate to each section and monitor staining closely.
13. As soon as the sections develop, immerse slides in dH<sub>2</sub>O.
14. Counterstain sections in hematoxylin.
15. Wash sections in dH<sub>2</sub>O two times for 5 minutes each.
16. Dehydrate sections.
17. Mount coverslips.