

NB100-56733 Protocol**Immunohistochemistry (Paraffin) Protocol**

Langerin/CD207 Antibody: https://www.novusbio.com/products/langerin-cd207-antibody_nb100-56733
 DeParaffinization

Materials Required

- Tissue array slide - 100%, 95%, and 75% ethanol
- Xylene

Method

1. Incubate in a dry oven at 62C for 1 hour. Slides should be maintained in a vertical orientation to allow complete removal of the paraffin.
2. Dewax slides in xylene for 5 x 4 minutes.
3. Hydrate slides in 100%, 95%, and 75% ethanol for 2 x 3 minutes each.
4. Immerse slides in tap water for 5 minutes.

Suggested Antigen Retrieval Protocol

The following procedure are a suggestion only. Other protocols can be used on the array slides.

Materials Required

- Tissue array slide - Phosphate buffered saline (pH 7.6)
- Citrate buffer (0.01 M, pH 6.0) - Microwave oven (700 W)

Method 1 (Microwave)

1. Immerse slides into citrate buffer (0.01 M, pH 6.0).
2. Microwave (700 W or high) for 5 min, add citrate buffer if necessary.
3. Microwave (medium) for 5 min, add citrate buffer if necessary.
4. Microwave (low) for 5 min.
5. Immerse in cold PBS.

Method 2 (Autoclave/Pressure Cooker)

1. Immerse slides in citrate buffer.
2. Incubate in a pressure cooker for 2 min at 95C.
3. Cool to room temperature.
4. Wash slides in PBS for 3 x 5 min.

Method 3 (Enzyme treatment)

1. Incubate slides with pronase [0.05% (w/v) in PBS] or trypsin [0.05% (v/v) in PBS] or pepsin [0.05% (v/v) in 2 N HCl] at 37oC (incubation time should be adjusted according to the antibody).
2. Wash slides in PBS for 3 x 5 min.

Method 4 (Hot bath)

1. Heat citrate buffer (1mM EDTA, pH8.0 or 0.01M sodium citrate buffer, pH6.0) to about 950C.
2. Place slides in the buffer for 10-15 min.

Immunostain**Materials Required**

1. Slides

2. Phosphate buffered saline (pH 7.6)
3. Hydrogen peroxide
4. Primary antibody
5. Blocking serum (normal serum)
6. Biotinylated secondary antibody
7. ABC reagent (6, 7, and 8 are included in Vectastain Elite ABC kit)
8. Diaminobenzidine
9. Meyer's hematoxylin
10. Permount

Method

1. Incubate in a dry oven at 62C for 1 hour. Slides should be maintained in a vertical orientation to allow complete removal of the paraffin.
2. Dewax slides in xylene for 5 x 4 minutes.
3. Hydrate slides in 100%, 95%, and 75% ethanol for 2 x 3 minutes each.
4. Immerse slides in tap water for 5 minutes
5. Antigen retrieval method (optional).
6. Quenching of endogenous peroxidase (optional)
 - a. immerse slides in 3% hydrogen peroxide solution for 6 minutes.
 - b. wash slides in PBS for 3 x 5 minutes.
7. Incubate slides with blocking serum (1:50) for 30 min.*
8. Blot excess serum from section, and incubate with primary Ab. Suggested incubation time may vary between antibodies:

mAb 2 hours at room temperature or overnight at 4C.

pAb: 1 ~ 1.5 hours at room temperature.
9. Wash slides in PBS for 3 x 5 minutes.
10. Incubate slides with biotin-conjugated secondary Ab for 30 min.*
11. Wash slides in PBS for 3 x 5 minutes.
12. Incubate slides with Avidin-Biotin Complexes for 30 min.*
13. Wash slides in PBS for 3 x 5 minutes.
14. Incubate slides in fresh DAB solution for 2 minutes. (We use DAB solution in Vector DAB/Ni substrate kit).**
15. Stop the reaction by washing in tap water.
16. Counterstain in Meyer's hematoxylin for 10 seconds.
17. Dehydrate slides in 75%, 80%, 95% and 100% ethanol
18. Clear slides in xylene 4 X 5 minutes.
19. Mount cover slide with Permount.

* Blocking serum, secondary antibody and avidin-biotin-peroxidase complexes are included in most of the immunostaining kit. Our lab uses the ABC kit from Vector Lab (Vectastain Elite ABC kit).

** We use DAB solution in Vector/DAB/Ni substrate kit (Vector Labs., Cat. SK-4100).