

NB300-122 Protocol

Immunohistochemistry protocol specific for TRPV1 Antibody (NB300-122)

Immunohistochemistry protocol specific for TRPV1 Antibody (NB300-122):

https://www.novusbio.com/products/vanilloid-r1-trpv1-antibody_nb300-122

Procedure Guide for NB 300-122 Polyclonal Anti-Vanilloid Receptor I

Immunohistochemistry Procedure

- 1) Animals were perfused transcardially with calcium-free Tyrode's solution, followed by fixative, and finally with 10% sucrose in PBS. Antiserum was used on perfusion-fixed tissues.
- 2) Tissues were dissected and stored in PBS/10% sucrose, overnight.
- 3) Frozen tissues were sectioned and mounted on gelatin-coated slides.
- 4) Slide-mounted tissue was processed for indirect immunofluorescence. Then the slides were incubated with blocking buffer for 1 hour at room temperature.
- 5) NB300-122 (anti-Vanilloid Receptor I) was diluted in blocking buffer to 1:2000.
- 6) Blocking buffer was removed and diluted NB300-122 was added. The primary antibody incubation was done at 4 degrees C for 18-24 hours.
- 7) Slides were rinsed 3 times and then incubated with appropriate secondary antibody for 1 hour at room temperature.
- 8) Slides were rinsed 3 times.
- 9) Slides were coverslipped.
- 10) Staining was examined by fluorescent microscopy.