



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NB300-201](http://www.novusbio.com/NB300-201)

## NB300-201 Protocol

# Serum protocol for Neurokinin B Antibody (NB300-201)

[[URL:[https://www.novusbio.com/products/neurokinin-b-antibody\\_nb300-201](https://www.novusbio.com/products/neurokinin-b-antibody_nb300-201)]] [[Caption:Neurokinin B Antibody]]  
Immunohistochemistry

1. Rats were perfused with 4% paraformaldehyde in 0.1M PBS.
2. The brains were then post-fixed in 4% paraformaldehyde and sucrose protected.
3. The treated sections were blocked in 0.1M Tris buffer containing 0.1% Triton X-100 and 0.25% BSA for 1 hour.
4. Anti-NKB (NB 300-201) was diluted in the blocking buffer and incubated with the sections overnight.
5. A goat anti-rabbit IgG conjugated to Alexa 488 was diluted in the blocking buffer and incubated with the sections overnight.

NOTE: Sections were made on a sledge microtome and stored in cryoprotectant solution at -20 degrees Celsius until ready to use.