



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/NB200-414](http://www.novusbio.com/NB200-414)

## NB200-414 Protocol

### Immunocytochemistry/Immunofluorescence Protocol for Influenza A H1N1 Nucleoprotein Antibody (NB200-414)

---

Immunocytochemistry/Immunofluorescence Protocol for Influenza A H1N1 Nucleoprotein antibody (NB200-414):

- 1) Cell monolayer was grown on a glass and infected with the virus for 2 days at 37C.
- 2) Glass was dried and the cells were fixed with cold acetone at 4C for 10 min.
- 3) NB 200-414 was introduced in concentrations from 1 to 100 ug/ml in PBS and incubated for 30 min in RT.
- 4) Glass was washed with PBS several times and incubated with goat anti-mouse IgG labeled with FITC.