



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB200-174

NB200-174 Protocol

Immunocytochemistry/Immunofluorescence Protocol for p19ARF Antibody (NB200-174)

[[URL:https://www.novusbio.com/products/p19arf-cdkn2a-antibody-5-c3-1_nb200-174]][[Caption:p19ARF Antibody]]
Immunofluorescent

1. Wild type mouse embryo fibroblasts (MEF) at passage 5 and NIH3T3 cells, which have deleted the Arf gene were fixed in methanol/acetone, and probed with anti-p19Arf (NB 200-174).
2. Cells were then probed with a fluorescently labeled anti-rat IgG secondary antibody.
3. Cells were also stained with Hoechst dye to reveal the position of nuclei.