



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/NB200-174](http://www.novusbio.com/NB200-174)

## NB200-174 Protocol

### Immunohistochemistry protocol for p19ARF/CDKN2A Antibody (NB200-174)

[[URL:[https://www.novusbio.com/products/p19arf-cdkn2a-antibody-5-c3-1\\_nb200-174](https://www.novusbio.com/products/p19arf-cdkn2a-antibody-5-c3-1_nb200-174)]][[Caption:p19ARF/CDKN2A Antibody]]

Immunohistochemistry

1. Sections from a wild type mouse testis taken at P11 were probed with anti-p19Arf (NB 200-174) [green], after antigen retrieval.
2. Sections were stained with DAPI to reveal the position of cell nuclei [blue].
3. p19Arf staining can be seen in subnuclear structures (presumably nucleoli) in cells at the periphery of the testicular tubules.