



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB100-428

NB100-428 Protocol

Immunohistochemistry protocol for FIH-1/HIF-1AN Antibody (NB100-428)

[[URL:https://www.novusbio.com/products/fih-1-hif-1an-antibody_nb100-428]][[Caption:FIH-1/HIF-1AN Antibody]]

1. Prepared 18 micron frozen rat brain tissue sections (post fixed with 10% formalin for 30 minutes).
2. Dilute the primary antibody 1:100 and incubated with diluted FIH antibody (cat# NB 100-428) at 4C, overnight.
3. Diluted the secondary antibody, biotinylated anti-rabbit IgG, according to the manufacturer's guidelines and incubated overnight at 4C.
4. Diluted the tertiary antibody, HRP-conjugated streptavidin, according to the manufacturer's guidelines and incubated overnight at 4C.
5. Visualized with DAB oxidation by HRP.