



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-00178

NBP1-00178 Protocol

Immunohistochemistry protocol for DIO2 Antibody (NBP1-00178)

Immunohistochemistry (NBP1-00178): https://www.novusbio.com/products/dio2-antibody_nbp1-00178

Specimen Preparation: Formalin-fixed, paraffin-embedded tissues are suitable for use with this primary antibody.

Deparaffinization: Deparaffinize slides using xylene or xylene alternative and graded alcohols.

Antibody Dilution: If using the concentrate format of this product, dilute the antibody 1:100. The dilutions are estimates; actual results may differ because of variability in methods and protocols.

Antigen Retrieval: Boil tissue section in EDTA buffer for 10 min followed by cooling at room temperature for 20 min.

Primary Antibody Incubation: Incubate for 10 minutes at room temperature.

Slide Washing: Slides must be washed in between steps. Rinse slides with PBS/0.05% Tween.

Visualization: Detect the antibody as instructed by the instructions provided with the visualization system.