



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-19048

NBP1-19048 Protocol

Immunohistochemistry protocol for Lubricin/PRG4 Antibody (NBP1-19048)

IHC Protocol (NBP1-19048): https://www.novusbio.com/products/lubricin-prg4-antibody_nbp1-19048
For use in IHC-Paraffin: Deparaffinize sections in xylene and graded ethanols. After deparaffinization, digest sections with Proteinase K for six minutes at room temperature followed by incubation in protein block serum free for ten minutes at room temperature.