



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-33779

NBP1-33779 Protocol

Western Blot protocol for p53 Antibody (NBP1-33779)

Western Blot (NBP1-33779): https://www.novusbio.com/products/p53-antibody-s11-k_nbp1-33779

- Wash buffer: 1x Tris Buffered Saline (TBS); 0.2% Tween-20
- Blocking buffer: 1xTBS; 0.2% Tween 20; 5% nonfat dry milk

For western blots, incubate the membrane with antibody diluted in blocking buffer 2 hours at room temperature.