



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-28570

NBP1-28570 Protocol

Western Blot protocol for Cytokeratin 19 Antibody (NBP1-28570)

NBP1-28570 Western blot Protocol (NBP1-28570): https://www.novusbio.com/products/cytokeratin-19-antibody-sb39g_nbp1-28570

Run protein on a 14% gel, using HCC38 & HT-29 cell lysate at 20 ug/well. (No sonication). Dilute NBP1-28570 to 2 ug/mL in 5% milk/TBST. 6mL/strip, incubated at RT on shaker for 1 hour. Dilute anti-mouse secondary antibody at 1:4K in 5% milk/TBST, 6mL/strip, incubate at RT on shaker for 1 hour.