



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB110-61646

NB110-61646 Protocol

Western Blot Protocol for Nucleophosmin Antibody (NB110-61646)

NPM1 Antibody: https://www.novusbio.com/products/npm1-antibody_nb110-61646
Western Blot

1. Run 50 ug of cytoplasmic fraction into each lane of an 12% SDS-PAGE.
2. Block the membrane.
3. Dilute the primary anti-mNPM1 (NB110-61646) in 1% NFDM and incubate for 2 hours at room temperature.
4. Wash the membrane.
5. Dilute the secondary anti-rabbit IgG, conjugated to HRP in 1% NFDM and incubate for 1 hour at room temperature.
6. Develop the blot.