



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB200-314

NB200-314 Protocol

Western Blot protocol for hnRNP M1-M4 Antibody (NB200-314)

hnRNP M1-M4 Antibody (1D8): https://www.novusbio.com/products/hnrnp-m1-m4-antibody-1d8_nb200-314
Western Blot Procedure Guide for NB 200-314 Monoclonal Anti-RNP M1-M4

- 1) Wet Nitrocellulose membrane with PBS + NP40 [PN].
- 2) Pour off PN.
- 3) Dilute anti-RNP M1-M4 (catalog# NB 200-314) in 5%NFDM + NP40.
- 4) Incubate the primary antibody with the membrane for 1 hour, at room temperature (RT), gently rocking.
- 5) Wash 1x with PN for 10 minutes. Wash 2x with PN for 5 minutes, each.
- 6) Dilute anti-mouse-HRP (Amersham) in 5%NFDM + NP40 at 1:5,000.
- 7) Incubate the secondary antibody with the membrane for 45 minutes, at RT, gently rocking.
- 8) Wash 1x with PN for 10 minutes. Wash 2x with PN for 5 minutes, each.
- 9) Wash 1x with PBS for 5 minutes.
- 10) Develop using Amersham ECL components.

****NOTE:** HeLa nuclear cell extracts can be used as a positive control for this antibody.