



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NB200-315](http://www.novusbio.com/NB200-315)

## NB200-315 Protocol

### Western Blot protocol specific for hnRNP M3-M4 Antibody (NB200-315)

hnRNP M Antibody (2A6): [https://www.novusbio.com/products/hnrnp-m-antibody-2a6\\_nb200-315](https://www.novusbio.com/products/hnrnp-m-antibody-2a6_nb200-315)  
Western Blot Procedure

- 1) Wet Nitrocellulose membrane with PBS + NP40 [PN].
- 2) Pour off PN.
- 3) Dilute anti-RNP M3/M4 (catalog# NB 200-315) 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.