

NB200-316 Protocol

Western Blot Protocol for CUGBP1 Antibody (NB200-316)

[[URL:https://www.novusbio.com/products/cugbp1-celf1-antibody-3b1_nb200-316]][[Caption:CUGBP1/CELF1 Antibody (3B1)]]

Procedure Guide for NB 200-316 Monoclonal Anti-CUG

Western Blot Procedure

- 1) Wet Nitrocellulose membrane with PBS + NP40 [PN].
- 2) Pour off PN.
- 3) Dilute anti-CUG (catalog# NB 200-316) 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.