

Results

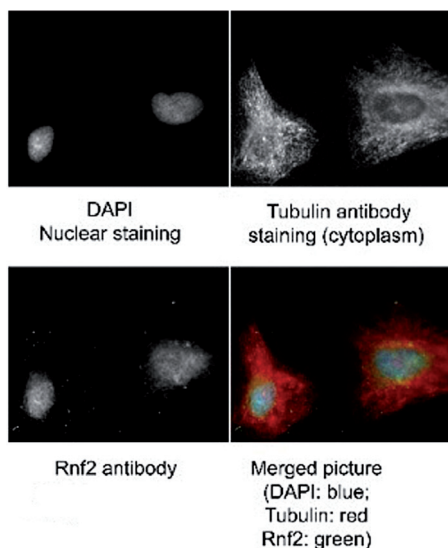


Figure 1. RING1B antibody Immunofluorescence results

Immunofluorescence of RING1B antibody. Tissue: human HeLa cells. Fixation in methanol and blocked with 0.2% fish scale gelatin for 1 hour at 25°C. Primary antibody at 1:300 for 20 minutes at 25°C. Secondary antibody: Alexa Fluor®488-conjugated Donkey anti-goat IgG secondary antibody at 1:500 for 45 min at RT. Localization: RING1B is nuclear and occasionally cytoplasmic. Staining: RING1B (RNF2) as green signal, Tubulin cytoplasm staining red, and DAPI (blue) nuclear counterstain.

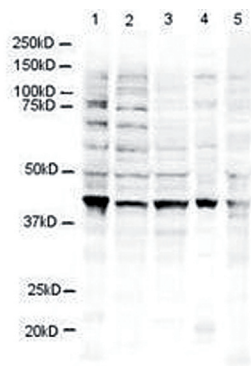


Figure 2. RING1B antibody western blot results

Western blot using the Diagenode RING1B antibody shows detection of a 38 kDa band corresponding to human RING1B in 3T3 (lane 1), U937 (lane 2), Jurkat (lane 3), mouse brain (lane 4) and CHO-K1 (lane 5) cell lysates. Approximately 20 µg of lysate was run on a SDS-PAGE and transferred onto nitrocellulose followed by reaction with a 1:500 dilution of RING1B antibody incubated at room temperature.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 10, 2014