

Results

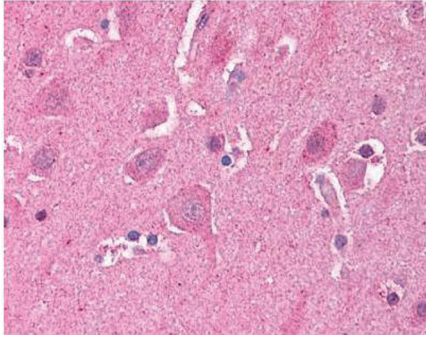


Figure 1. LDB2 antibody immunohistochemistry results

The Diagenode LDB2 antibody was used at 5 µg/ml to detect LDB2 in human brain cortex tissue. The image shows the localization of antibody as the precipitated red signal, with a hematoxylin purple nuclear counter stain. Tissue was formalin-fixed and paraffin embedded.

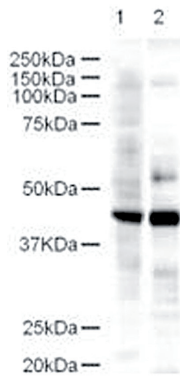


Figure 2. LDB2 antibody Western blot results

Western blot using the Diagenode LDB2 antibody shows detection of a 43-kDa band corresponding to LDB2 in a lysates prepared from human kidney (lane 1) and mouse spleen (lane 2) tissues. Approximately 18 µg of lysate was run on a SDS-PAGE and transferred onto nitrocellulose followed by reaction with a 1:500 dilution of LDB2 antibody.

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