

LDB2 polyclonal antibody - Classic

Cat. No. C15410271

Type: Polyclonal

Source: Rabbit

Lot #: 001

Size: 100 µg

Concentration: 1.06 µg/µl

Specificity: Mouse, human, chimpanzee, dog, rat, chicken

Purity: Affinity purified

Storage: Store at -20°C; for long storage, store at -80°C.
Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Applications

	Suggested dilution	Results
Immunohistochemistry	1:200	Figure 1
Western blot	1:500 - 1:2,000	Figure 2

Target description

It is suggested that LIM domain binding proteins (LDB1, LDB2 and LDB3) act synergistically to enhance transcriptional efficiency by acting as co-factors for LIM homeodomain and Otx class transcription factors both of which have essential roles in development. LDB2 is a close homologue of LDB1. LDBs are homologs of the Drosophila Chip protein.

Results

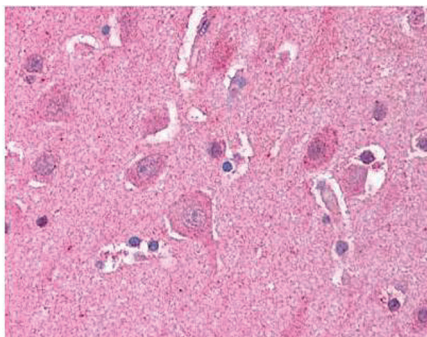


Figure 1. LDB2 antibody immunohistochemistry results

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.

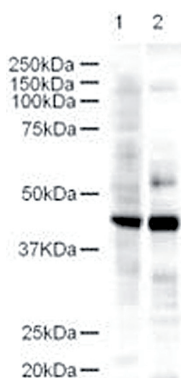


Figure 2. LDB2 antibody Western blot results

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