

LDB1 polyclonal antibody - Classic

Cat. No. C15410269

Type: Polyclonal

Source: Rabbit

Lot #: 001

Size: 100 µg

Concentration: 1.80 µg/µl

Specificity: Human, mouse, rat

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
ELISA	1:425,000	
Western blot	1:500 - 1:3,000	Figure 1

Target description

LDB1 is also known as CLIM 2, LIM Domain Binding 1, NLI and Nuclear LIM Domain Interactor. The LIM-domain binding protein binds to the LIM domain of LIM homeodomain proteins which are transcriptional regulators of development. Nuclear LIM interactor (NLI) / LIM domain-binding protein 1 (LDB1) is located in the nuclei of neuronal cells during development, it is co-expressed with Isl1 in early motor neuron differentiation and has a suggested role in the Isl1 dependent development of motor neurons. It is suggested that these proteins 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.

Results

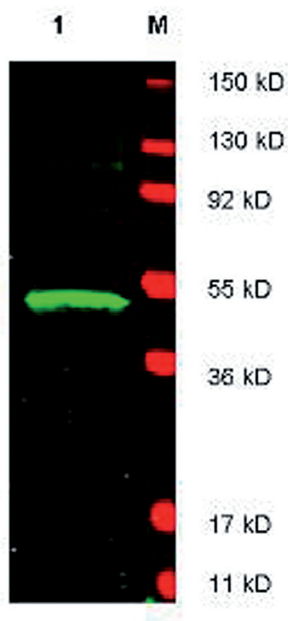


Figure 1. LDB1 antibody Western blot results

Western blot using the Diagenode LDB1 antibody shows detection of LDB1 protein in Jurkat whole cell lysate. Approximately 30 μ g of lysate was loaded prior to separation and transfer to nitrocellulose. Primary antibody was used at a 1:1,800 dilution. Molecular weight estimation was made by comparison to prestained MW markers in (lane M)

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