

## IGF2BP2 antibody (other name: VICKZ2, IMP-2, IMP2)

**Cat. No.** C15410350

**Type:** Polyclonal

**Source:** Rabbit

**Lot:** 001

**Size:** 100 µg

**Concentration:** 1 µg/µl

**Specificity:** Mouse: positive.

Other species: not tested.

**Purity:** Affinity purified polyclonal antibody.

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

**Storage buffer:** PBS containing 50% glycerol, does not contain a preservative.

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

**Description:** Polyclonal antibody raised in rabbit against human IGF2BP2 (Insulin Like Growth Factor 2 mRNA Binding Protein 2), using KLH-conjugated synthetic peptide from the C-terminal part of the protein.

### Applications

| Applications     | Suggested dilution                 | References |
|------------------|------------------------------------|------------|
| RIP              | 15 µg per 10 <sup>7</sup> cells    | Fig 1      |
| Western Blotting | 1:1,000                            | Fig 2      |
| IP               | 5 µg per 2.5x10 <sup>6</sup> cells | Fig 3      |

### Target description

IGF2BP2 (UniProtKB/Swiss-Prot entry Q9Y6M1) is a member of the insulin-like growth factor 2 mRNA-binding protein family. It binds to the 5' UTR of the mRNA of certain genes, IGF2, ACTB and MYC, thereby regulating translation. Binding of IGF2BP2 recruits the target mRNA to cytoplasmic protein-RNA complexes (mRNPs) which allows for mRNA transport and transient storage. Further, it modulates the rate of translation and protects the mRNA from degradation. IGF2BP2 plays an important role in metabolism and has been associated with susceptibility to diabetes.

#### Diagenode sa. BELGIUM | EUROPE

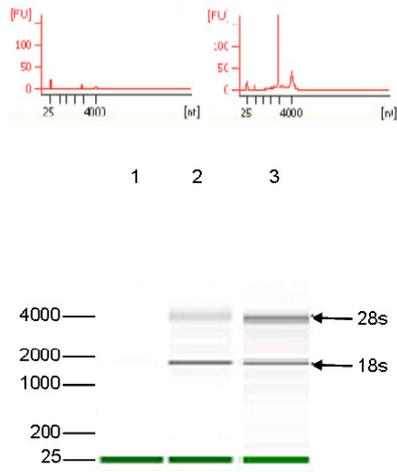
LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

#### 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

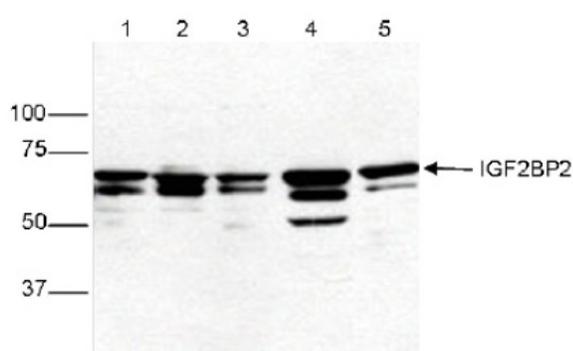
Last update: March, 2018

## Results



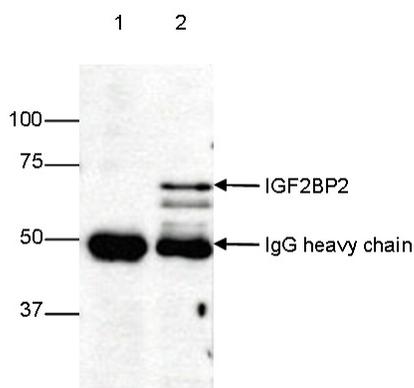
**Figure 1. Immunoprecipitation using the Diagenode antibody directed against IGF2BP2**

Immunoprecipitation was performed on total RNA isolated from 10 million K562 cells using 15 µg of the Diagenode antibody against IGF2BP2 (cat. No. C15410350) or with an equal amount of rabbit IgG, used as a negative control. The immunoprecipitated RNA was subsequently analysed on a Bioanalyzer. Figure 1 shows the Bioanalyzer profile obtained with the negative control (upper left) and the IGF2BP2 antibody (upper right). The lower figure shows the gel image for the negative IgG control, the IGF2BP2 antibody and the input (lane 1, 2 and 3 respectively). The marker (in bp) is shown on the left, the position of the 28s and 18s ribosomal RNA is indicated on the right.



**Figure 2. Western blot analysis using the Diagenode antibody directed against IGF2BP2**

Whole cell extracts from K562, 293T, HeLa, NIH3T3 and Rat1 cells (lanes 1, 2, 3, 4 and 5, respectively) were analysed by Western blot using the Diagenode antibody against IGF2BP2 (cat. No. C15410350) diluted 1:1,000 in TBS containing 1% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 3. Immunoprecipitation using the Diagenode antibody directed against IGF2BP2**

Immunoprecipitation was performed on whole cell extracts from K562 cells using 5 µg of the Diagenode monoclonal antibody against IGF2BP2 (cat. No. C15410350, lane 2). An equal amount of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated IGF2BP2 protein was subsequently detected by western blot with the IGF2BP2 antibody as described above.