

## IGF2BP1 monoclonal antibody

**Cat. No. C15200237**

Type: Monoclonal	Specificity: Human, mouse: positive. Other species: not tested.
Size: 100 µg	Isotype: IgG2a
Concentration: 1 µg/µl	Host: Mouse
Lot No.: 001	Purity: Protein A purified monoclonal antibody.
Storage buffer: PBS containing 50% glycerol, does not contain a preservative.	Storage conditions: Store at -20°C.
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Last Data Sheet Update: February 23, 2018

### Description

**Other names:** CRD-BP, VICKZ1, IMP-1, CRDBP, ZBP-1, IMP1, ZBP1

Monoclonal antibody raised in mouse against human IGF2BP1 (Insulin Like Growth Factor 2 mRNA Binding Protein 1), using the full length recombinant protein.

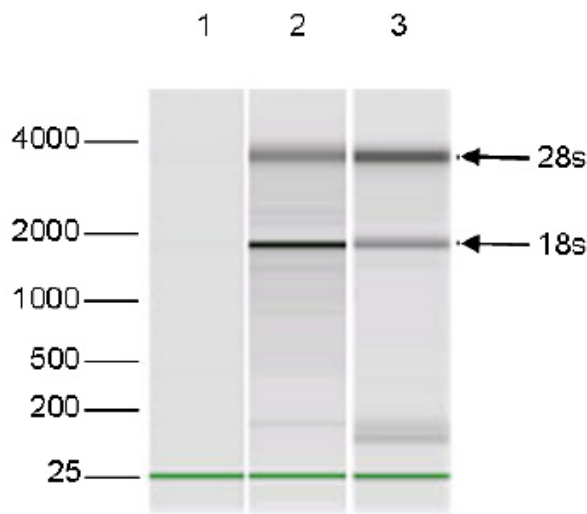
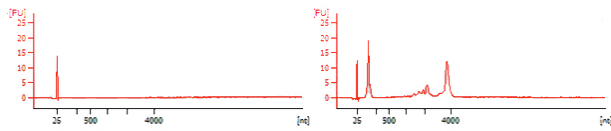
### Applications

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 5x10 <sup>6</sup> cells	Fig 3

### Target Description

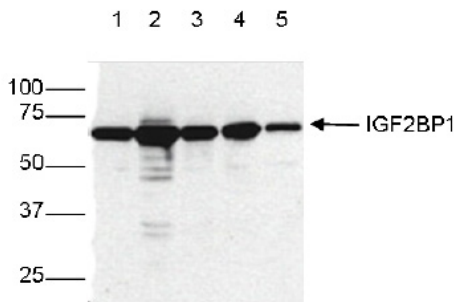
IGF2BP1 (UniProtKB/Swiss-Prot entry Q9NZI8) is a member of the insulin-like growth factor 2 mRNA-binding protein family. It binds to the mRNAs of certain genes, including insulin-like growth factor 2, beta-actin and beta-transducin repeat-containing protein, thereby regulating translation.

**Validation data**



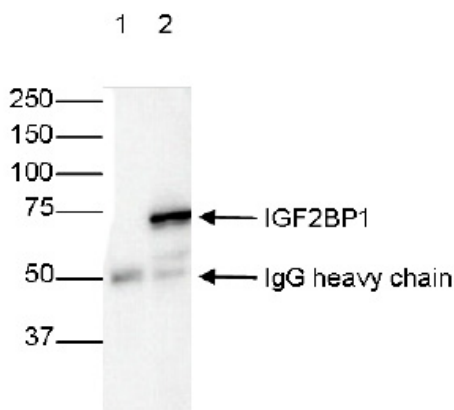
**Figure 1. Immunoprecipitation using the Diagenode monoclonal antibody directed against IGF2BP1**

Immunoprecipitation was performed on total RNA isolated from 10 million HeLa cells using 15 µg of the Diagenode antibody against IGF2BP1 (cat. No. C15200237) or with an equal amount of mouse IgG2a, 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 IGF2BP1 antibody (upper right). The lower figure shows the gel image for the negative IgG2a control, the IGF2BP1 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 monoclonal antibody directed against IGF2BP1**

Whole cell extracts from HeLa, 293T, HL-60, NIH3T3 and MEF cells (lanes 1, 2, 3, 4 and 5, respectively) were analysed by Western blot using the Diagenode monoclonal antibody against IGF2BP1 (cat. No. C15200237) 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 monoclonal antibody directed against IGF2BP1**

Immunoprecipitation was performed on whole cell extracts from K562 cells using 5 µg of the Diagenode monoclonal antibody against IGF2BP1 (cat. No. C15200237, lane 2). An equal amount of mouse IgG2a was used as a negative control (lane 1). The immunoprecipitated IGF2BP1 protein was subsequently detected by western blot with the IGF2BP1 antibody as described above.