

## IGF2BP3 polyclonal antibody

Cat. No. C15410352

Type: Polyclonal	Specificity: Human, mouse: positive. Other species: not tested.
Size: 100 µg	Isotype: NA
Concentration: 1 µg/µl	Source: Rabbit
Lot No.: 001	Purity: Affinity purified polyclonal 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: March 27, 2018

### Description

**Other names:** VICKZ3, IMP-3, IMP3, KOC, KOC1, HKOC, CT98

Polyclonal antibody raised in rabbit against human IGF2BP3 (Insulin Like Growth Factor 2 mRNA Binding Protein 3), using a KLH-conjugated synthetic peptide from the C-terminal part of the protein.

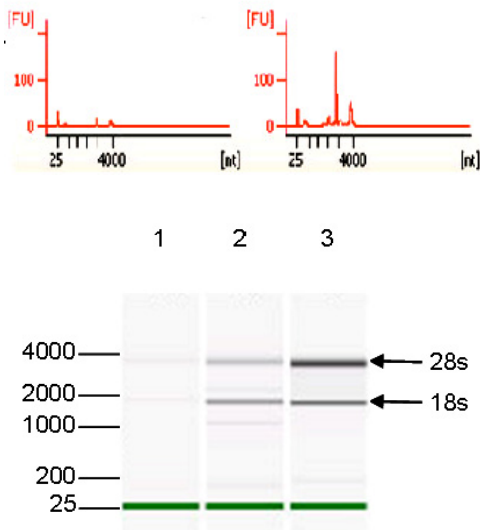
### Applications

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

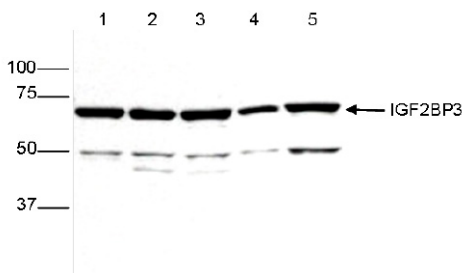
IGF2BP3 (UniProtKB/Swiss-Prot entry O00425) is a member of the insulin-like growth factor 2 mRNA-binding protein family. It binds to the 5' UTR of the IGF2 and some other mRNA. It may repress translation of IGF2 during late development. Binding of IGF2BP3 recruits the target mRNA to cytoplasmic protein-RNA complexes (mRNPs) which allows for mRNA transport and transient storage. Further, it modulates the rate of translation and protects the mRNA from degradation. IGF2BP also stabilizes the 3' UTR of CD44 mRNA, promoting cell adhesion and invadopodia formation in cancer cells.

**Validation data**



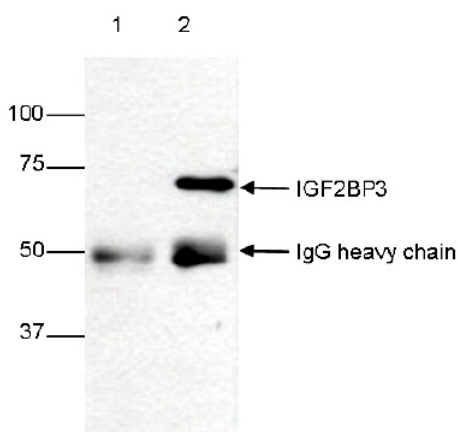
**Figure 1. Immunoprecipitation using the Diagenode antibody directed against IGF2BP3**

Immunoprecipitation was performed on total RNA isolated from 10 million K562 cells using 15 µg of the Diagenode antibody against IGF2BP3 (Cat. No. C15410352) 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 IGF2BP3 antibody (upper right). The lower figure shows the gel image for the negative IgG control, the IGF2BP3 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 IGF2BP3**

Whole cell extracts from K562, 293T, HeLa, Jurkat and NIH3T3 cells (lanes 1, 2, 3, 4 and 5, respectively) were analysed by Western blot using the Diagenode antibody against IGF2BP3 (Cat. No. C15410352) diluted 1:1,000 in PBS 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 IGF2BP3**

Immunoprecipitation was performed on whole cell extracts from K562 cells using 5 µg of the Diagenode monoclonal antibody against IGF2BP3 (Cat. No. C15410352, lane 2). An equal amount of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated IGF2BP3 protein was subsequently detected by western blot with the IGF2BP3 antibody as described above.