

## ELAVL1 monoclonal antibody

Cat. No. C15200238

Type: Monoclonal	Specificity: Human, mouse, rat, hamster: positive. Other species: not tested.
Size: 100 µg	Isotype: IgG2a
Concentration: 1 µg/µl	Source: Mouse
Lot No.: 001	Purity: 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: March 28, 2018

### Description

**Other names:** ELAV, ELAV1, MeIG, HUR, Hua

Monoclonal antibody raised in mouse against human ELAVL1 (ELAV Like RNA Binding Protein 1).

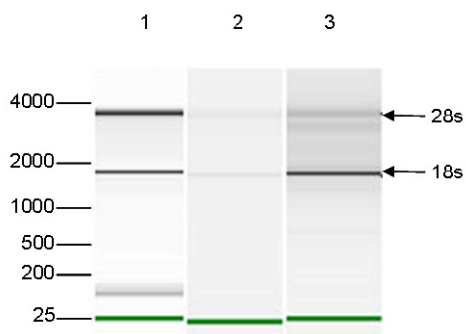
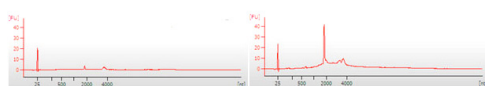
### Applications

Applications	Suggested dilution	References
RIP	15 µg per 10 <sup>7</sup> cells	Fig 1
Western blotting	1:1,000	Fig 2
IP	1.5 µg per 1.5x10 <sup>5</sup> cells	Fig 3

### Target Description

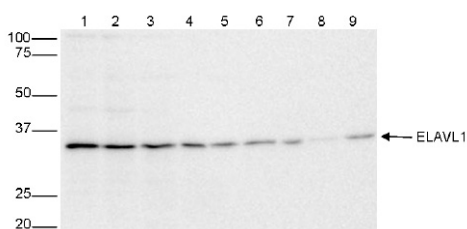
ELAVL1 (UniProtKB/Swiss-Prot entry Q15717) is a member of the ELAVL RNA-binding protein family. It selectively binds to poly-U elements and AU-rich elements (AREs) in the 3-UTR of target mRNAs, thereby stabilizing the mRNA. ELAVL1 has been implicated in a variety of biological processes including embryonic stem cells (ESCs) differentiation and has been linked to a number of diseases, including cancer. It is highly expressed in many cancers, and could be potentially useful in cancer diagnosis, prognosis, and therapy.

**Validation data**



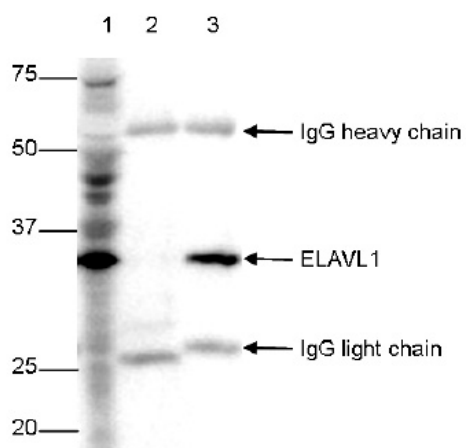
**Figure 1. Immunoprecipitation using the Diagenode monoclonal antibody directed against ELAVL1**

Immunoprecipitation was performed on total RNA isolated from 10 million HeLa cells using 15 µg of the Diagenode monoclonal antibody against ELAVL1 (Cat. No. C15200238) 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 ELAVL1 antibody (upper right). The lower figure shows the gel image for the input, the negative IgG control and the ELAVL1 antibody (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 ELAVL1**

Whole cell extracts from 293T, HeLa, K562, Jurkat, NIH3T3, WR19L, Rat1, PC12 and CHO cells (lanes 1 to 9, respectively) were analysed by Western blot using the Diagenode monoclonal antibody against ELAVL1 (Cat. No. C15200238) 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 monoclonal antibody directed against ELAVL1**

Immunoprecipitation was performed on whole cell extracts from HeLa cells using 1.5 µg of the Diagenode monoclonal antibody against ELAVL1 (Cat. No. C15200238, lane 3). An equal amount of mouse IgG2a was used as a negative control (lane 2). Lane 1 shows the input. The immunoprecipitated ELAVL1 protein was subsequently detected by western blot with the ELAVL1 antibody as described above.