

## ELAVL3 polyclonal antibody

**Cat. No. C15410348**

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: April 9, 2018

### Description

**Other names:** HUC, HUCL, PLE21

Polyclonal antibody raised in rabbit against human ELAVL3 (ELAV Like RNA Binding Protein 3), using a KLH-conjugated synthetic peptide from the N-terminal part of the protein.

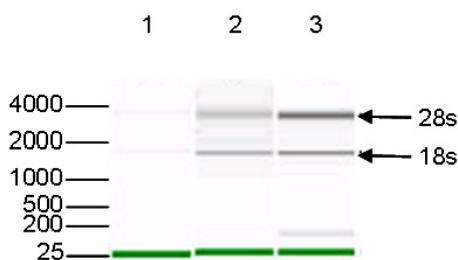
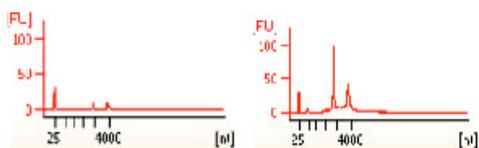
### Applications

Applications	Suggested dilution	References
RIP	15 µg per 6x10 <sup>6</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

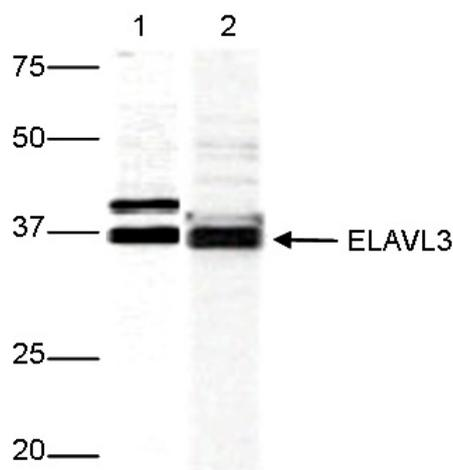
ELAVL3 (UniProtKB/Swiss-Prot entry Q14576) is a neural-specific member of the ELAVL RNA-binding protein family. It selectively binds to AU-rich sequences in several mRNA's including VEGF mRNA. The observation that ELAVL3 is recognized by the anti-Hu antibody present in sera from patients with paraneoplastic encephalomyelitis and sensory neuronopathy suggests it is involved in neuronal differentiation and maintenance.

**Validation data**



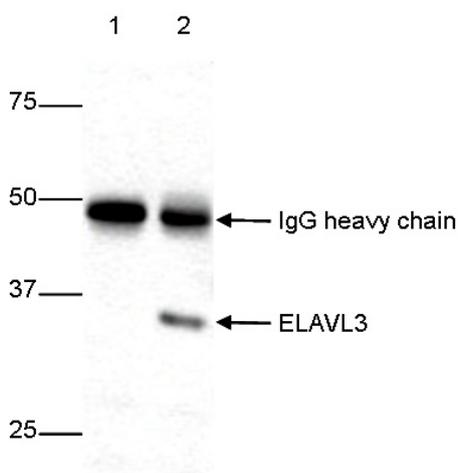
**Figure 1. Immunoprecipitation using the Diagenode antibody directed against ELAVL3**

Immunoprecipitation was performed on total RNA isolated from 6 million SK-N-SH cells using 15 µg of the Diagenode antibody against ELAVL3 (Cat. No. C15410348) 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 ELAVL3 antibody (upper right). The lower figure shows the gel image for the negative IgG control, the ELAVL3 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 ELAVL3**

Whole cell extracts from SK-N-SH cells (lane 1) and from P19 cells treated with retinoic acid (lane 2) were analysed by Western blot using the Diagenode antibody against ELAVL3 (Cat. No. C15410348) 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 ELAVL3**

Immunoprecipitation was performed on whole cell extracts from SK-N-SH using 5 µg of the Diagenode antibody against ELAVL3 (Cat. No. C15410348, lane 2). An equal amount of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated ELAVL3 protein was subsequently detected by western blot with the ELAVL3 antibody as described above.