

PABPC4 polyclonal antibody

Cat. No. C15410358

Type: Polyclonal	Specificity: Human, mouse, rat: 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: May 22, 2018

Description

Other names: PABP4, PABP-4, APP1, APP-1, IPABP

Polyclonal antibody raised in rabbit against human PABPC4 (Poly(A) Binding Protein Cytoplasmic 4), using a KLH-conjugated synthetic peptide from the central region of the protein.

Applications

Applications	Suggested dilution	References
RIP	15 µg per 10 ⁷ cells	Fig 1
Western blotting	1:1,000	Fig 2
IP	5 µg per 5x10 ⁶ cells	Fig 3

Target Description

PABPC4 (UniProtKB/Swiss-Prot entry Q13310) is a Poly(A)-binding protein. These proteins bind to the poly(A) tail present at the 3-prime ends of most eukaryotic mRNAs and may be involved in cytoplasmic regulatory processes of mRNA metabolism. PABPC4 probably also binds to cytoplasmic RNA sequences other than poly(A) in vivo. Further, it was identified as an antigen (APP1, activated-platelet protein-1), expressed on thrombin-activated platelets and may be involved in the regulation of protein translation in platelets and megakaryocytes.

Validation data

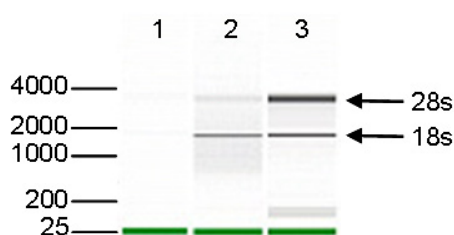
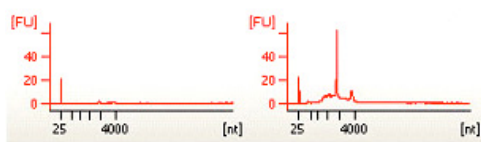


Figure 1. Immunoprecipitation using the Diagenode antibody directed against PABPC4

Immunoprecipitation was performed on total RNA isolated from 10 million 293T cells using 15 µg of the Diagenode antibody against PABPC4 (Cat. No. C15410358) 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 PABPC4 antibody (upper right). The lower figure shows the gel image for the negative IgG control, the PABPC4 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.

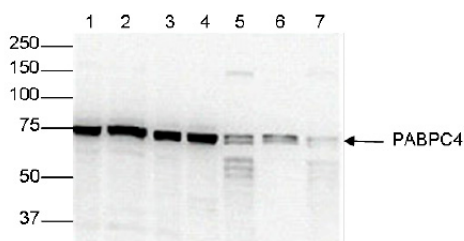


Figure 2. Western blot analysis using the Diagenode antibody directed against PABPC4

Whole cell extracts from 293T, HeLa, K562, Jurkat, NIH3T3, WR19L and Rat1 cells (lanes 1 to 7, respectively) were analysed by Western blot using the Diagenode antibody against PABPC4 (Cat. No. C15410358) 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.

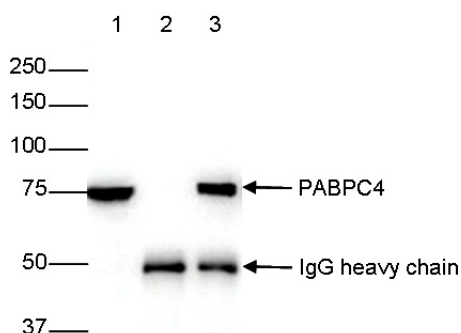


Figure 3. Immunoprecipitation using the Diagenode antibody directed against PABPC4

Immunoprecipitation was performed on whole cell extracts from 293T cells using 5 µg of the Diagenode antibody against PABPC4 (Cat. No. C15410358, lane 3). An equal amount of rabbit IgG was used as a negative control (lane 2). The immunoprecipitated PABPC4 protein was subsequently detected by western blot with the PABPC4 antibody as described above. The input is shown in lane 1.