

PRODUCT NAME		
PAPD4 polyclonal antibody		
Other names: GLD2		
Cat. No. C15310130 (CS-130-100)	Type: Polyclonal	Size: 100 µl
Lot #: A400-001	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human PAPD4 (PAP associated domain containing 4), using a KLH-conjugated synthetic peptide containing a sequence from the N-terminal part of the protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:1,000	Fig 1
Western blotting	1:1,000	Fig 2

Purity: Whole antiserum from rabbit containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Last data sheet update: March 18, 2010

Target description

PAPD4 (UniProtKB/Swiss-Prot entry Q6PIY7) is an RNA polymerase that generates the poly(A) tail of specific RNAs by adding AMP monomers to the 3'-end. In contrast to the canonical nuclear poly(A) RNA polymerase, this activity is restricted to selected cytoplasmic mRNAs.

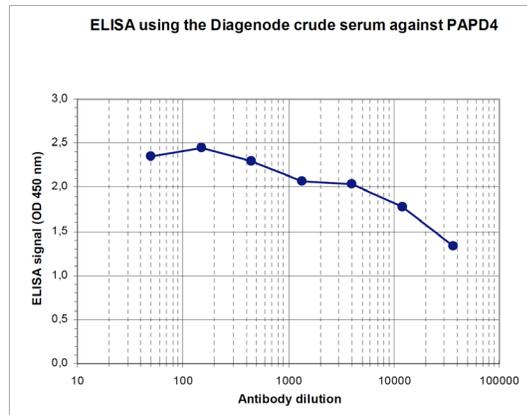


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human PAPD4 (Cat. No. CS-130-100). The plates were coated with the peptide used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:107,000.

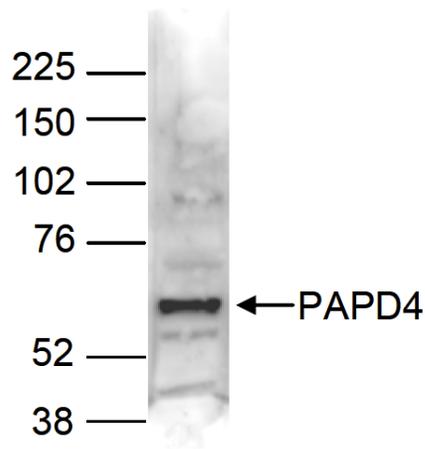


Figure 2
Western blot analysis using the Diagenode antibody directed against PAPD4

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against PAPD4 (Cat. No. CS-130-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right (expected size: 56 kDa); the marker (in kDa) is shown on the left.