

PRODUCT NAME		
PAPD5 polyclonal antibody		
Other names: TRF4-2		
Cat. No. C15310134 (CS-134-100)	Type: Polyclonal	Size: 100 µl
Lot #: A322-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human PAPD5 (PAP associated domain containing 5), using a KLH-conjugated synthetic peptide containing a sequence from the C-terminal part of the protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1

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

PAPD5 (UniProtKB/Swiss-Prot entry Q8NDF8) is an RNA polymerase which is involved in DNA repair. It also seems to have a function in sister chromatid cohesion.

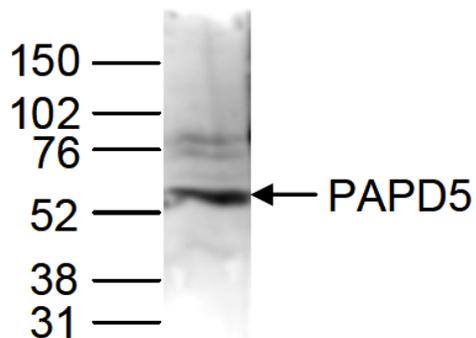


Figure 1

Western blot analysis using the Diagenode antibody directed against PAPD5

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against PAPD5 (Cat. No. CS-134-100) 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: 63 kDa); the marker (in kDa) is shown on the left.