

PRODUCT NAME		
Med26 polyclonal antibody		
Other names: Transcriptional coactivator CRSP70, ARC70		
Cat. No. <b>C15410073</b> (pAb-073-050)	Type: Polyclonal	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

**Description:** Polyclonal antibody raised in rabbit against human Med26 (Mediator of RNA polymerase II transcription subunit 26), 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
Western blotting	1:1,000	Fig 1

**Purity:** Protein G purified, polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

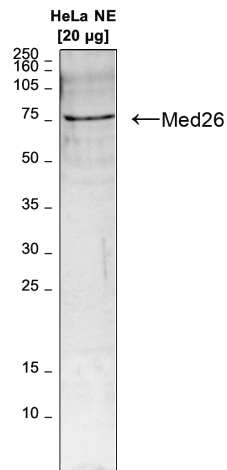
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

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

**Last data sheet update:** March 17, 2010

**Target description**

Med26 [UniProt/Swiss-Prot entry O95402] is a component of the Mediator complex, a coactivator involved in the regulation of transcription of nearly all RNA polymerase II-dependent genes. The Mediator complex forms a bridge between gene-specific regulatory proteins and the RNA polymerase II transcription machinery. It is recruited to promoters by direct interactions with these regulatory proteins and serves as a scaffold for the assembly of a functional pre-initiation complex with RNA polymerase II and general transcription factors.



**Figure 1**

**Western blot analysis using the Diagenode antibody directed against Med26**

Western blot was performed on nuclear extracts from HeLa cells (20  $\mu$ g) with the Diagenode antibody against human Med26 (Cat. No. pAb-073-050), diluted 1:1,000 in TBS-Tween containing 5% skimmed milk (Figure 2). The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.