

PRODUCT NAME		
CCDC101 polyclonal antibody		
Other names: SGF29		
Cat. No. C15310122 (CS-122-100)	Type: Polyclonal	Size: 100 µl
Lot #: A509-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human CCDC101 (coiled-coil domain containing 101), using two KLH-conjugated synthetic peptides containing a sequence from the central and from the C-terminal part of the protein, respectively.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:200	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 23, 2010

Target description

CCDC101 (UniProt/Swiss-Prot entry Q96ES7) associates with the histone acetyl transferase complexes TFIIIC-HAT and STAGA and may therefore be involved in transcriptional regulation. It has been shown to activate c-Myc target gene expression and may play a role in c-Myc-mediated tumorigenesis.

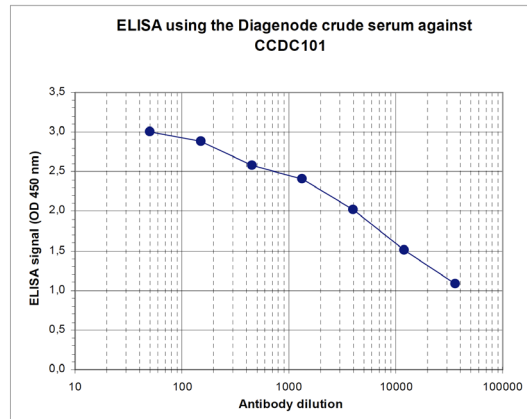


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human CCDC101 [Cat. No. CS-122-100]. The plates were coated with the peptides 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 1:13,700.

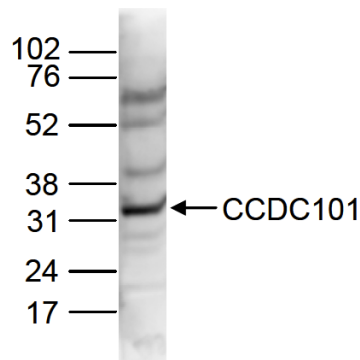


Figure 2
Western blot analysis using the Diagenode antibody directed against CCDC101

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against CCDC101 [Cat. No. CS-122-100] diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest (expected size: 33 kDa) is indicated on the right; the marker (in kDa) is shown on the left.