

PRODUCT NAME		
Wdr5 polyclonal antibody		
Other names: BIG3, SWD3		
<b>Cat. No.</b> C15310101 (CS-101-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A266-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse Wdr5 (WD (tryptophan-aspartate) repeat domain protein 5), using two KLH-conjugated synthetic peptides containing an amino acid sequence from the central part of the protein (1).

**Specificity:** Mouse: positive  
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:1,000	Fig 2, (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.

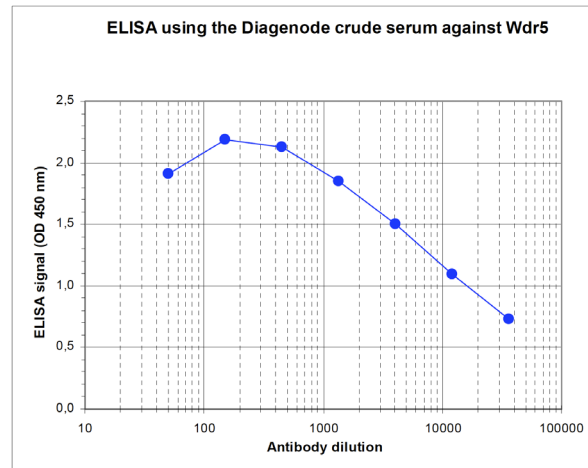
**References:**

(1) Peptide design by Andrea Kranz, Western blot analysis by Heike Petzold and Andrea Kranz, BIOTEC, Dept. of Genomics, Prof. F. Stewart, TU Dresden, Tatzberg 47/49, 01307 Dresden, Germany

**Last data sheet update:** April 7, 2010

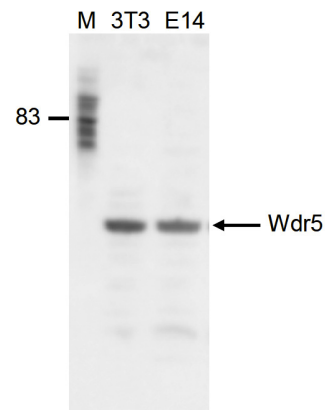
**Target description**

Wdr5 (UniProtKB/Swiss-Prot entry P61964) belongs to the family of WD repeat proteins which are involved in different cellular processes such as cell cycle progression, signal transduction, apoptosis, and gene regulation. It is a component of the Set1A and Set1B histone H3 methylation complexes. These complexes methylate lysine 4 of H3, thereby activating gene transcription. Wdr5 interacts with H3 dimethyl K4, but not with tri or mono methylated H3K4.



**Figure 1**  
**Determination of the titer**

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse Wdr5 [Cat. No. CS-101-100]. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:12,300.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against Wdr5 (1)**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Wdr5 [Cat. No. CS-101-100], diluted 1:1,000 in BSA/PBS-Tween. The molecular weight marker (M, in kDa) is shown on the left; the location of the protein of interest (36 kDa) is indicated on the right.