

## RBBP5 polyclonal antibody

**Other names:** RBBP-5, RBQ3, SWD1

**Cat. No.** C15410342

**Type:** Polyclonal **ChIP grade/ChIP-seq grade**

**Source:** Rabbit

**Lot #:** A300-109A2

**Size:** 100 µl

**Concentration:** 1 µg/µl

**Specificity:** Human: positive

Other species: not tested

**Purity:** Affinity purified polyclonal antibody in Tris-citrate buffer containing 0.09% azide.

**Storage:** Store at 4°C.

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

**Description:** Polyclonal antibody raised in rabbit against human RBBP5 (RB Binding Protein 5), using a synthetic peptide containing a sequence from the C-terminus of the protein<sup>1</sup>.

### Applications

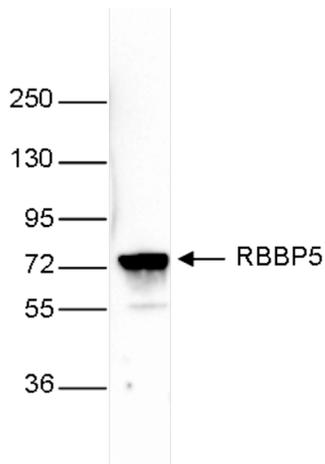
Applications	Suggested dilution	References
ChIP*	2 µg per ChIP	Fig 1, 2
Western blotting	1:1,000	Fig 3
IP	10 µg per IP	Fig 4
IHC	1:1,000	Fig 5

\*Please note that the optimal antibody amount per ChIP should be determined by the end-user. We recommend testing 1-5 µg per ChIP.

### Target description

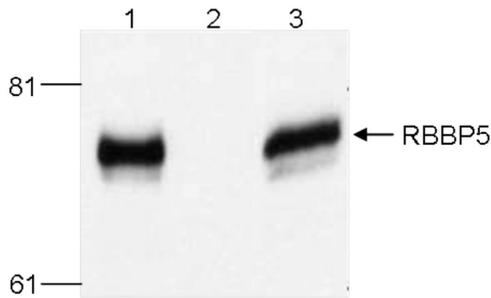
RBBP5 (UniProtKB/Swiss-Prot entry Q15291) belongs to the highly conserved family of WD-repeat proteins. It's part of the MLL1/MLL complex and is as such involved in mono-, di- and trimethylation at lysine 4 of histone H3, a specific tag for epigenetic transcriptional activation. RBBP5 binds to the retinoblastoma protein thereby playing a role in the regulation of cell proliferation. It also plays a crucial role in the differentiation potential in embryonic stem (ES) cells.





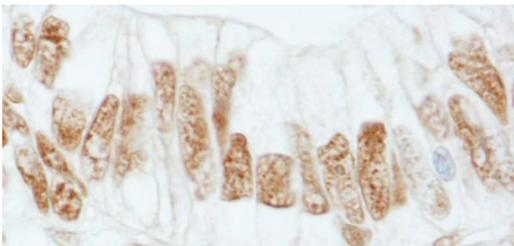
**Figure 3. Western blot analysis using the Diagenode antibody directed against RBBP5**

Whole cell extracts from HeLa cells were analysed by Western blot using the Diagenode antibody against RBBP5 (Cat. No. C15410342) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 4. Immunoprecipitation using the Diagenode antibody directed against RBBP5**

Immunoprecipitation was performed on whole cell extracts from 293T cells using 10 µg of the Diagenode antibody against RBBP5 (Cat. No. C15410342, lane 1). An equal amount of rabbit IgG was used as a negative control (lane 2). The immunoprecipitated CBX2 protein was detected by western blot with the RBBP5 antibody diluted 1:500. Lane 3 shows the input (15% of the IP).



**Figure 5. Immunohistochemistry using the Diagenode antibody directed against RBBP5**

FFPE sections of human colon carcinoma tissue were analysed by Immunohistochemistry with the Diagenode antibody against RBBP5 (Cat. No. C15410342, lane 1) diluted 1:1,000.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** May 25, 2017