

## Smc3K112/113ac (yeast) polyclonal antibody

**Other names:** BAM, CSPG6, SMC3L1, HCAP, CDLS3

**Cat. No.** C15310229 (CS-PA009-100)

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** A1012-001

**Size:** 100 µl

**Concentration:** not determined

**Specificity:** Yeast: positive / Other species: not tested

**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.

**Description:** Polyclonal antibody raised in rabbit against yeast Smc3 (structural maintenance of chromosomes 3) acetylated at lysines 112 and 113, using a KLH-conjugated synthetic peptide.

### Applications

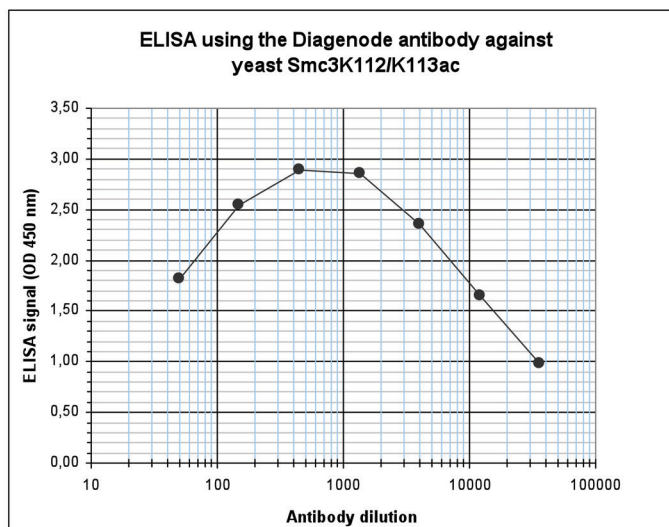
	Suggested dilution	Results
ELISA	1:1,000	Fig 1

\*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:1,000 dilution

### Target description

SMC3 (UniProtKB/Swiss-Prot entry Q9UQE7) is a central component of the cohesion complex, which is required for chromosome cohesion during the cell cycle. The cohesion complex may form a large proteinaceous ring within which sister chromatids can be trapped. At anaphase, the complex is cleaved and dissociates from chromatin, allowing sister chromatids to segregate. Cohesion is coupled to DNA replication and is involved in DNA repair. The cohesin complex plays also an important role in spindle pole assembly during mitosis and in chromosome movement. Acetylation of SMC3 is required for the establishment of cohesion during cell division.

## Results



**Figure 1. Determination of the antibody titer**

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against yeast Smc3K112/113ac (cat. No. CS-PA009-100). The plates were coated with the peptide 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 1:16,900.

### **Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - 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:** September 4, 2012