

<b>PRODUCT NAME</b> G9a polyclonal antibody		
Other names: EHMT2, BAT8, KMT1C, NG36		
<b>Cat. No.</b> C15310096 (CS-096-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A262-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse G9a (Protein G9a), 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	1:500 – 1:1,000	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.

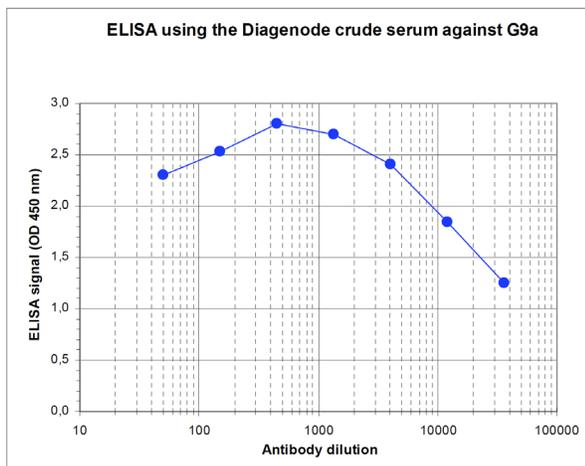
**References:**

(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

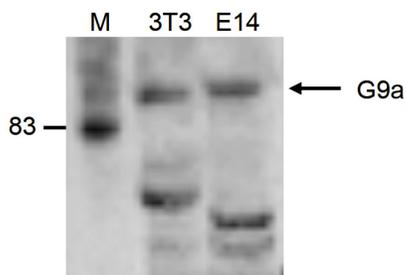
**Target description**

G9a [UniProtKB/Swiss-Prot entry Q96KQ7] is a member of the Set 1 family of histone methyltransferases. It preferentially methylates K9 and K27 of histone H3. Methylation of H3K9 represents a specific tag for epigenetic transcriptional repression by recruiting HP1 proteins to methylated histones. G9a is a part of the E2F6.com-1 complex.



**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 G9a [Cat. No. CS-096-100]. The wells were coated with the peptides used for immunisation of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:30,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against G9a (1)**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) cells with the Diagenode antibody against mouse G9a [Cat. No. CS-096-100], 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 (138 kDa) is indicated on the right.