

PRODUCT NAME		
FKBP51 polyclonal antibody		
Other names: FKBP5, FKBP54, AIG6, PPlase, Rotamase		
<b>Cat. No.</b> C15310110 (CS-110-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A371-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse FKBP51 [FK506 Binding Protein 51], 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	1:500	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.

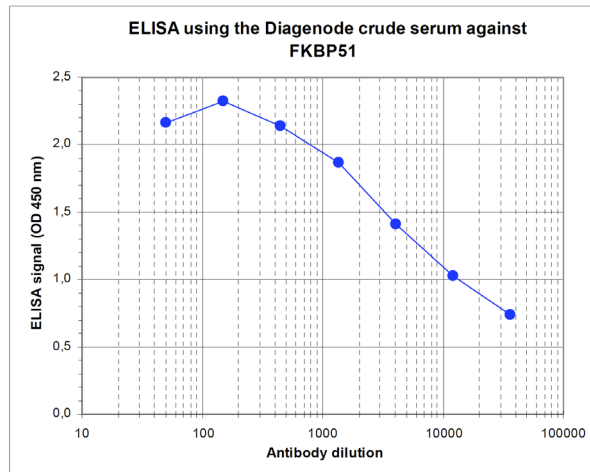
**References:**

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

**Last data sheet update:** April 6, 2010

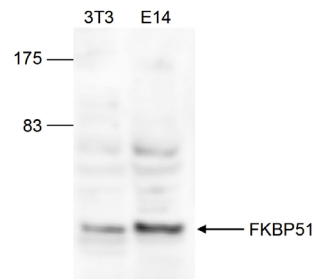
**Target description**

FKBP51 [UniProtKB/Swiss-Prot entry Q13451] is a member of the immunophilin protein family, which plays a role in immunoregulation and basic cellular processes involving protein folding and trafficking. It acts as a receptor for FK506 and rapamycin, two immunosuppressants which inhibit T-cell proliferation. FKBP51 is thought to mediate calcineurin inhibition and interacts with progesterone receptor complexes.



**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 FKBP51 (Cat. No. CS-110-100). The plates were coated with a mix of 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 1:9,400.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against mouse FKBP51 (1)**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse FKBP51 (Cat. No. CS-110-100), diluted 1:500 in BSA/PBS-Tween. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest (expected size: 51 kDa) is indicated on the right.