

PRODUCT NAME		
SPI1 polyclonal antibody		
Other names: PU.1, SFPI1, OF, SPI-A, SPI-1		
<b>Cat. No.</b> C15310120 (CS-120-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A516-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse SPI1 (spleen focus forming virus [SFFV] proviral integration oncogene), using two KLH-conjugated synthetic peptides containing a sequence from the N-terminus and from the C-terminus of the protein, respectively.

**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

**Purity:** Whole antiserum from rabbit containing 0.05% azide.

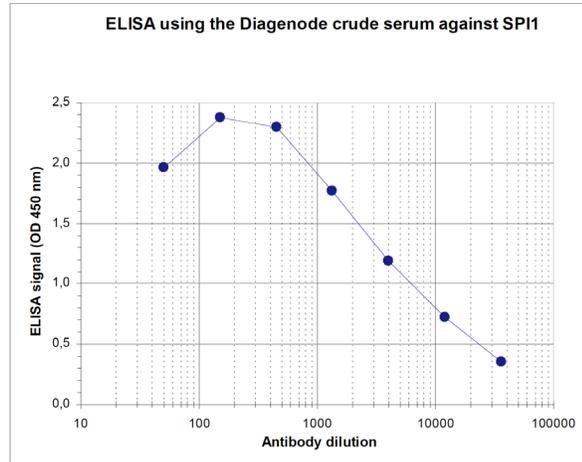
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

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

**Last data sheet update:** March 19, 2010

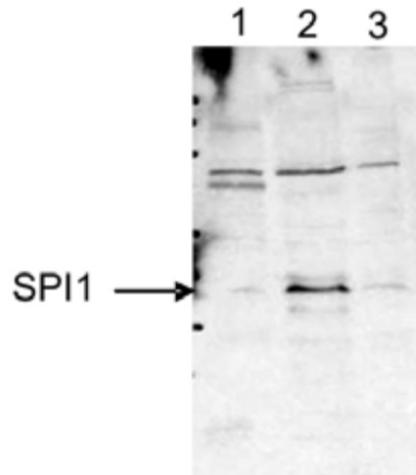
**Target description**

SPI1 (UniProt/Swiss-Prot entry P17947) is an ETS-domain transcription factor that binds to a purine-rich DNA sequence (5'-GAGGAA-3') found near the promoters of target genes, also known as the PU-box. SPI1 is a transcriptional activator that may be involved in myeloid and B-lymphoid cell development. Further, SPI1 may also be important for the regulation of splicing of target genes.



**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 SPI1 (Cat. No. CS-120-100). The plates were coated with 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 1:2,600.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against SPI1**

Western blot was performed using and the Diagenode antibody against mouse SPI1 (Cat. No. CS-120-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. Figure 2 shows the results for mouse fibroblasts (3T3, lane 1), macrophages (RAW, lane 2) and pre-B cells (HAFTL, lane 3). The position of the protein of interest (expected size: 31 kDa) is indicated on the left. Fibroblasts do not express SPI1, whereas macrophages show a high and pre-B cells a low expression of SPI1.

Western blot performed by Maribel Parra, Center for Genomic Regulation, Barcelona, Spain.