

Flp polyclonal antibody

Cat. No. C15310169

Type: Polyclonal	Specificity:
Size: 100 µl	Isotype: NA
Concentration: not determined	Host: Rabbit
Lot No.: A280-004	Purity: Whole antiserum from rabbit containing 0.05% azide.
Storage buffer: NA	Storage conditions: 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: February 12, 2021

Description

Polyclonal antibody raised in rabbit against Flp recombinase using 3 KLH-conjugated synthetic peptides located at the N-terminal part of the protein ⁽¹⁾.

Applications

Applications	Suggested dilution	References
ELISA	1:500	Fig 1
Western Blotting	1:500	Fig 2, (1)

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

Target Description

Flp recombinase originates from *Saccharomyces cerevisiae* and recognizes FRT sites in the genome. It catalyzes recombination between two FRT sites, thereby excising the intervening DNA sequence. Flp recombinase is used as a tool for the generation of transgenic animals.

Validation data

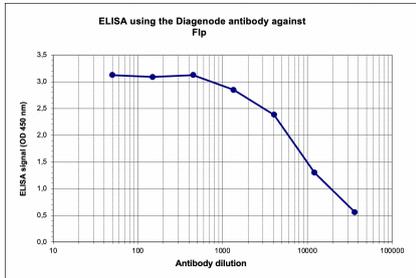


Figure 1. Determination of the antibody titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against Flp (cat. No. C15310169) in antigen coated wells. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:9,000.

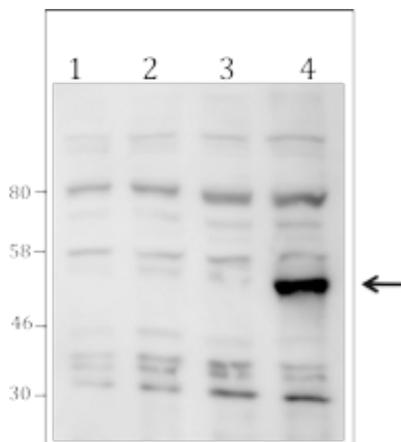


Figure 2. Western blot analysis using the Diagenode antibody directed against Flp⁽¹⁾

Western blot was performed on whole cell lysates from untransfected 293 cells (lane 1), or 293 cells transfected with Cre (lane 2), Dre (lane 3) or Flp (lane 4) with the Diagenode antibody against Flp (cat. No. C15310169), diluted 1:500 in BSA/PBS-Tween. The molecular weight marker (in KD) is shown on the left; the location of the protein of interest (expected size: 48 KD) is indicated on the right.