



6405 Mira Mesa Blvd, Ste 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

## Data Sheet

### Anti-Cre polyclonal antibody Catalog #: 25215

<b>Lot #:</b> 220819	<b>Host Species:</b> Rabbit
<b>Conc.:</b> N/A	<b>Species Reactivity:</b> NA
<b>Size:</b> 100 µl	<b>Immunogen:</b> Synthetic peptide
<b>Clonality:</b> Polyclonal	<b>Purification:</b> Whole serum

**Description:** Polyclonal antibody raised in rabbit against Cre recombinase using a KLH-conjugated synthetic peptide located at the C-terminal part of the protein.

**Background:** Cre recombinase originates from bacteriophage P1 and recognizes loxP sites in the genome. It catalyzes recombination between two loxP sites, thereby excising the intervening DNA sequence. Cre recombinase is used as a tool for the generation of transgenic animals

**Formulation:** Whole antiserum from rabbit containing 0.05% azide

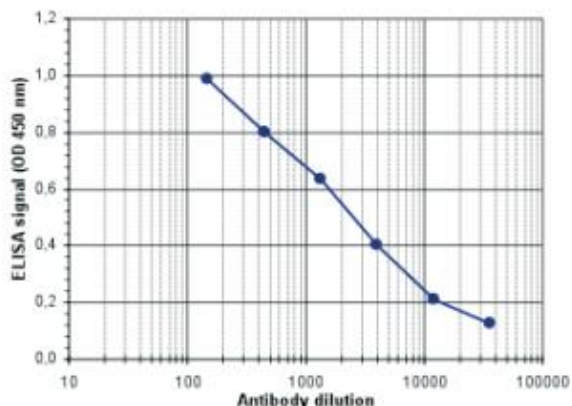
**Applications:** ELISA (1:100), WB (1:500)

**Storage/Stability:** Store at -80°C for up to 2 years. Centrifuge after first thaw to maximize product recovery. Aliquot to avoid repeated freeze/thaw cycles. Aliquots may be stored at -20°C for at least one month.

**Warnings:** Avoid freeze/thaw cycles

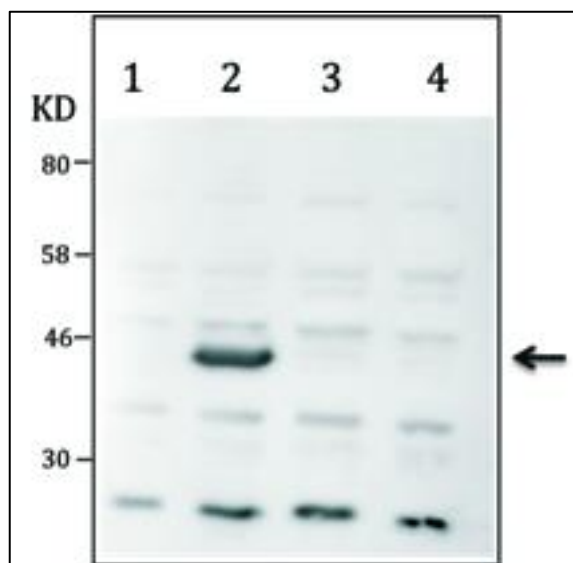
**Notes:** The optimal antibody concentration should be determined by the end-user.

**Quality Assurance:**



**Determination of the antibody titer.**

To determine the titer, an ELISA was performed using a serial dilution of the antibody directed against Cre (cat. No. 25215) in antigen coated wells. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:2,900.



**Western blot analysis using the antibody directed against Cre (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 antibody against Cre (cat. No. 25215), 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: 38 kDa) is indicated on the right.