

## PrP<sup>C</sup>His

**Genotyp: ARR (Codons 136,154,171), scrapie resistant ovine (sheep), recombinant, *E. coli***

Cat. No.	Amount
PR-376	50 µg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### **Avoid freeze / thaw cycles**

### **Form**

Supplied as lyophilized powder.

### **Purity**

> 95% by SDS-PAGE.

### **Description**

N-terminal His-tagged full-length mature PrP<sup>C</sup>. The protein may be reconstituted in detergent-containing buffers e.g. TX-100 (0.5%) or under mild denaturing conditions (e.g. 1.5 M guanidine-HCl or urea).

The His-PrP<sup>C</sup> is an ideal positive control in immunochemical detection of prions. It may be used directly for the production of antibodies in animals.

### **Selected References:**

- Wopfner et al. (1999) Analysis of 27 mammalian and 9 avian PrPs reveals high conservation of flexible regions of the prion protein. *J. Mol. Biol.* **289**:1163.
- Houston et al. (2003) Prion diseases: BSE in sheep bred for resistance to infection. *Nature.* **423**:498.
- Thorgeirsdottir et al. (1999) PrP gene polymorphism and natural scrapie in Icelandic sheep. *J. Gen. Virol.* **80**:2527.