

PrP^CHis

**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

Lyophilized. Lyophilized from 10mM Tris-HCl pH6.7.

Purity

> 95% by SDS-PAGE

Description

N-terminal His-tagged full-length mature PrP^C. 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^C 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.