

## PrP<sup>c</sup> His

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

Cat. No.	Amount
PR-377	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.  
The protein may be reconstituted indetergent-containing buffers e.g. 0.5% TX-100 or under mild denaturing conditions (e.g. 1.5 M guanidine-HCl or urea).

#### Purity

> 95% by SDS-PAGE

#### Description

N-terminal His-tagged full-length mature PrP<sup>c</sup>.  
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. His-PrP<sup>c</sup> (VRQ-Genotyp) is also a substrate of choice for studies of conversion of PrP<sup>c</sup> into PrP<sup>Res</sup> compared to the ARR Genotyp.

#### 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.