

**native PrP<sup>c</sup> (full length, residues 23-231)<sup>His</sup>**  
**(Prion Protein, cellular form)**  
**Human, Recombinant, *E. coli***

Cat. No.	Amount
PR-908	5 µg

Supplied as frozen liquid (50 µM PrP) in 10 mM sodium acetate, pH 4.5 and 0.02 % sodium azide.

N-terminal His-tagged full-length prion protein (PrP<sup>c</sup>).

The PrP<sup>c</sup>-like conformation of the protein after re-folding procedure was confirmed by high resolution NMR spectroscopy.

The native His tagged PrP<sup>c</sup> is the ideal tool for *in vitro/in vivo* investigations of the behaviour of cellular prions. It may be used for protein-protein interaction studies as well as for *in vitro/in vivo* conversion studies.

AVOID FREEZE/THAW CYCLES!

**For *in vitro* use only!**

**Purity:** >95% by SDS-PAGE

**Store:** -20 °C or -80 °C

## Selected References:

Prusiner S.B. (1998). Prions. *Proc. Natl. Acad. Sci. USA* **95**:13363.

Pan *et al.* (1993) Conversion of  $\alpha$ -helices into  $\beta$ -sheets features in the formation of the scrapie prion proteins. *Proc. Natl. Acad. Sci. USA* **90**:10962.

Lee *et al.* (1998) Complete genomic sequence and analysis of the prion protein gene region from three mammalian species. *Genome Res.* **8**:1022.

Bergstrom *et al.* (2005) Amidation and Structure Relaxation Abolish the Neurotoxicity of the Prion Peptide PrP106-126 in Vivo and in Vitro. *J. Biol. Chem.* **280**:23114.