

PrP^c (full length, residues 23-231)^{His}

Prion Protein, cellular form
human, recombinant, *E. coli*

Cat. No.	Amount
PR-902	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.

Molecular weight

25165 Da.

Purity

>90% by SDS-PAGE.

Description

Recombinant N-terminal His-tagged full-length prion protein. 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:

- Prusiner S.B. (1998) Prions. *Proc. Natl. Acad. Sci. USA* **95**:13363.
Pan et al. (1993) Conversion of α -helices into β -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.
Nixon R.R. (2005) Prion-associated Increases in Src-family Kinases. *J. Biol. Chem.* **280**:2455.
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.