

PrP^c (globular domain, residues 90-231)^{His}

Prion Protein, cellular form

human, recombinant, *E. coli*

| Cat. No. | Amount |
|----------|--------|
| PR-905 | 50 µg |

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

Avoid freeze / thaw cycles

Form

Lyophilized. Dialyzed before against 0.1 µM HCL.

Molecular Weight

18.5 kDa

Purity

>90% by SDS-PAGE

Description

N-terminal His-tagged prion protein globular domain (amino acid residues 90-231). The protein may be re-constituted 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-tagged 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 *et al.* (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.