

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

Prion Protein, cellular form

hamster, recombinant, *E. coli*

Cat. No.	Amount
PR-909	5 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 10 mM sodium acetate pH 4.5 and 0.02 % sodium azide.

Purity

>95% by SDS-PAGE

Description

N-terminal His-tagged full-length prion protein (PrP^c). The PrP^c-like conformation of the protein after refolding procedure was confirmed by high resolution NMR spectroscopy.

The native His tagged PrP^c 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.

Selected References:

- Prusiner (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.