

## PrP<sup>c</sup> (globular domain, residues 90-230)<sup>His</sup>

(Prion Protein, cellular form)

Mouse, Recombinant, *E. coli*

Cat. No.	Amount
PR-907	50 µg

Supplied as lyophilized powder.

N-terminal His-tagged prion protein globular domain (amino acid residues 90-230). 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-tagged 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.

AVOID FREEZE/THAW CYCLES!

**For in vitro use only!**

**Purity:** >90% by SDS-PAGE

**Store:** -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.