

PrP^c (full length, residues 23-231)^{His}
(Prion Protein, cellular form)
Hamster, Recombinant, *E. coli*

Cat. No.	Amount
PR-903	50 µg

Supplied as lyophilized powder.

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.

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 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.