

PrP^c (F209S)^{His}

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

Cat. No.	Amount
PR-372	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.

Purity

>95% by SDS-PAGE.

Description

N-terminal His-tagged mutant (F209S) mature PrP^c. 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 bovine PrP^c (F209S) is a homologue to the disease-associated human (F198S) point mutation. It may be used to investigate the effect of this mutation for protein stability in *in vitro*/*in vivo* conversion studies.

Selected References:

- Glockshuber R. (2001). Folding dynamics and energetics of recombinant prion proteins. *Adv. Prot. Chem.* **57**:83.
- 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.