

PrP^c (A144V)^{His}

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

Cat. No.	Amount
PR-374	50 µg

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

Avoid freeze / thaw cycles

Form

Lyophilized. Lyophilized from 10mM Tris-HCl pH6.7.

Purity

>95% by SDS-PAGE

Description

N-terminal His-tagged mutant (A144V) 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 (A144V) is a homologue to the scrapie-sensitive bovine mutant A136V. It may be used to investigate the effect of this mutation on protein stability in *in vitro*/*in vivo* conversion studies.

Selected References:

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

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.