

bovPrP^c cDNA [A144V] cDNA from bovine Prion Protein, cellular form

Cat. No.	Amount
PRD-374	10 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Supplied in 10 mM Tris-HCl, pH 8.0.

Description

BovPrPc cDNA wild type, encoding the full-length mature protein (amino acids 25-242).

ATG-start codon and different restriction sites for sub-cloning on request.

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.