

## Prion Protein Monoclonal Antibody - Pri 308



### ORDERING INFORMATION

**Catalogue Number:** A03200

**Size:** 200 µg

**Stability:** ≥2 years

**Storage:** -20°C

**Shipping:** Dry Ice

- **Host species:** Mouse
- **Antigen species:** This anti-prion Protein Monoclonal Antibody was raised against a synthetic peptide representative of the human 106-126 sequence,
- **Label:** Labelling upon request
- **Form:** lyophilized IgG
- **Clone:** Pri 308
- **Isotype:** IgG
- **Application:** EIA, WB
- **Description:**

Prion Protein (PrP) and namely its abnormal isoform, partially resistant to proteinase K (PrPres), is the only specific molecular marker of the Transmissible Spongiform Encephalopathies (TSEs) such as Bovine Spongiform Encephalopathie (BSE) or its human form, the New Variant of Creutzfeld-Jakob disease.

**For research laboratory  
use only.**

**Not for human  
diagnostic use.**

***Should you wish further information, do not hesitate to contact us.***

