

## Prion Protein Monoclonal Antibody - SAF 83



### ORDERING INFORMATION

**Catalogue Number:** A03207

**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 preparation of SAF (scrapie associated fibrils) from infected hamster brain.
- **Label:** Labelling upon request
- **Form:** lyophilized IgG
- **Clone:** SAF 83
- **Isotype:** IgG
- **Application:** EIA, FC, 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.***

