

Prion Protein Monoclonal Antibody - 8G8



ORDERING INFORMATION

Catalogue Number: A03220

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 the human recombinant protein.
- **Label:** Labelling upon request
- **Form:** lyophilized IgG
- **Clone:** 8G8
- **Isotype:** IgG
- **Application:** IHC
- **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.

