

Prion Protein Monoclonal Antibody - Sha31



ORDERING INFORMATION

Catalogue Number: A03213

Size: 200 µg

Storage: -20°C

Shipping: Dry Ice

- **CAS Number:** 0
- **Host species:** Mouse
- **Antigen species:** Protein Kinase treated and non-denaturated SAF from Syrian Hamster
- **Form:** Ascitic liquid lyophilized
- **Clone:** Sha 31
- **Isotype:** IgG1
- **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.
- **Articles:**
 - ▣ Krasemann S, Neumann M, Geissen M, Bodemer W, Kaup F-J, et al. (2010) Preclinical Deposition of Pathological Prion Protein in Muscle of Experimentally Infected Primates. PLoS ONE 5(11): e13906. doi:10.1371/journal.pone.0013906
 - ▣ Vulin J, Biacabe AG, Cazeau G, Calavas D, and Baron T Molecular typing of protease-resistant prion protein in transmissible spongiform encephalopathies of small ruminants, France, 2002-2009. Emerg Infect Dis, Jan 2011; 17(1): 55-63.

For research laboratory use only.

Not for human diagnostic use.

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

