

Product Information



Prion Protein Monoclonal Antibody - BAR 233

Cat No: A03223 - 200 µg

General Data

Shipping:	Dry Ice
Stability:	Store at -20°C; shelf life 2 years maximum after production
Formulation:	lyophilized IgG with BSA
Host:	Mouse
Antigen:	This anti-Prion Protein Monoclonal Antibody was raised against an ovine recombinant protein.
Clone:	BAR233
Isotype:	IgG2b
Label:	Labelling upon request



Application(s): Reconstitute the content of the vial in 1 mL of water.

For western blot analysis of PrP_C, dilute to a final concentration of 1 µg/mL and use BSA as blocking reagent. For EIA and flow cytometry, the optimal working must be determined empirically (currently between 0.1 and 1 µg/mL).

Specificity: PrP_C
(NA*) Hamster, (+) Mouse, (NA*) Bovine, (+) Ovine, (+) Human.

Special advice: GMO product: require notification of the end-consumers GMO Registration Number and the name of the person at the facility who is responsible for GMO Supervision & Safety

Product Overview

Prion Protein (PrP) and namely its abnormal isoform, partially resistant to proteinase K (PrP^{Sc}), 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. This antibody recognises the protein sequence within amino acids 141-152 (human numbering).

Scientific Literature

Féraudet C., et al. Screening of 145 Anti-PrP Monoclonal Antibodies for Their Capacity to Inhibit PrP^{Sc} Replication in Infected Cells. The journal of Biological Chemistry. Vol. 280. No. 12, March 25, pp. 11247-11258, 2005.

FP/17/16

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Pharma does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Pharma be liable for incidental, consequential or punitive damages.

Contact Bertin Pharma

10 bis avenue Ampère
F-78180 Montigny-le-Bretonneux
Tel.: +33 (0)139 306 036
bioreagent@bertinpharma.com
bioreagent.bertinpharma.com

Product Information



Prion Protein Monoclonal Antibody - BAR 233

Cat No: A03223 - 200 µg

Selective and Efficient Immunoprecipitation of the Disease-associated Form of the Prion Protein Can Be Mediated by Nonspecific Interactions between Monoclonal Antibodies and Scrapie-associated Fibrils. The Journal of Biological Chemistry. Vol. 279. No. 29, July 16, pp. 30143-30149, 2004.

FP/17/16

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Pharma does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Pharma be liable for incidental, consequential or punitive damages.

Contact Bertin Pharma

10 bis avenue Ampère
F-78180 Montigny-le-Bretonneux
Tel.: +33 (0)139 306 036
bioreagent@bertinpharma.com
bioreagent.bertinpharma.com