



## Prion Protein Monoclonal Antibody - SAF 61

Cat No: A03205 - 200 µg

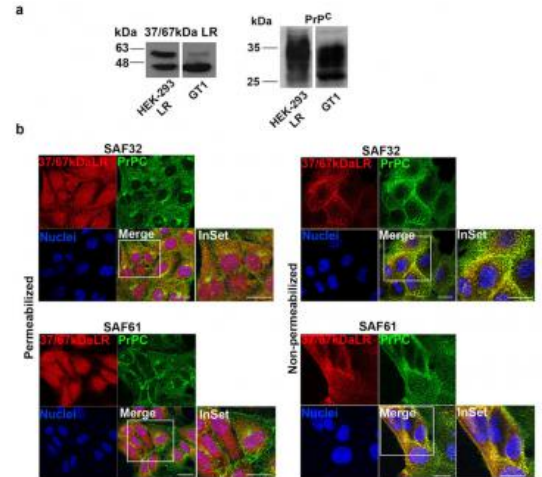
### General Data

<b>Shipping:</b>	Dry Ice
<b>Stability:</b>	Store at -20degrees;shelf life 2 years maximum after production
<b>Formulation:</b>	lyophilized IgG with BSA
<b>Host:</b>	Mouse
<b>Antigen:</b>	This anti-Prion Protein Monoclonal Antibody was raised against a preparation of SAF (scrapie associated fibrils) from infected hamster brain.
<b>Clone:</b>	SAF 61
<b>Isotype:</b>	IgG2a
<b>Label:</b>	Labelling upon request

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

For EIA and flow cytometry, the optimal working dilution must be determined empirically (currently between 0.1 and 1 µg/mL).

**Specificity:** PrPc  
(+) Hamster, (+) Mouse, (+) Bovine, (+) Ovine, (+) Human



### Product Overview

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.

This antibody recognises the protein sequence within amino acids 142-160 (human numbering).

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

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac  
10 bis avenue Ampère  
78180 Montigny le Bretonneux - France

Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>

<https://www.bertin-bioreagent.com>