

# Product Information



## SMRT<sup>Å</sup> Monoclonal Antibody

Cat No: G01038 - 100 ÅµL

### General Data

<b>Shipping:</b>	wet ice
<b>Stability:</b>	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
<b>Formulation:</b>	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
<b>Host:</b>	Mouse
<b>Antigen:</b>	recombinant human SMRT protein
<b>Clone:</b>	1000000
<b>Isotype:</b>	IgG2a, Å
<b>Application(s):</b>	ELISA Western Blot Immunofluorescence Recommended dilutions: 1/500-1/5000
<b>Specificity:</b>	human SMRT protein. Epitope between aa 982 and 1695



FP/19/20

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'Activit s du Pas du Lac  
10 bis avenue Amp re  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036

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