

## Data Sheet

### Anti-GST Tag

Mouse Monoclonal Antibody  
Anti-GST Tag (GST.B6 / G2R)  
Catalog #: 25009

#### Specificity:

Recognizes native and denatured forms of purified GST or GST fusion proteins

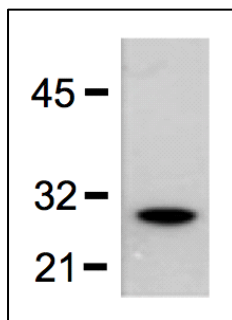
#### Immunogen:

Recombinant Glutathione-S-transferase protein (*S. japonicum*)

#### Applications:

Dot Blot, ELISA, ICC, IP, WB

#### Image:



**Fig.** 1:5000 (0.2µg/mL) Ab dilution probed against BL-21 bacteria lysate expressing the GST protein

#### Clonality | Host Species | Isotype:

GST.B6 / G2R | Mouse | IgG2a

#### Formulation:

1mg/mL in 10mM PBS, pH 7.2, with 0.05% NaN<sub>3</sub>

#### Sensitivity:

WB (with ECL): 10-50ng of purified GST or GST fusion proteins

#### Working Conditions:

WB (with ECL) 1:1000-3000 dilution (incubate for 1 hour at room temperature)  
ICC: 1:500-2000 dilution  
For best results with other assays (e.g.: Dot Blot, ELISA, IP, etc.), please determine optimal working dilution by titration test

#### Purification:

Protein A affinity chromatography from mouse ascites fluid

#### Storage and Stability:

Store at -20°C for up to 2 years. Centrifuge after first thaw to maximize product recovery. Aliquot to avoid repeated freeze-thaw cycles. Store aliquots at 4°C for 1-2 weeks

OUR PRODUCTS ARE FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)