

Anti-GST

Mouse Monoclonal Antibody

Catalog # **G52-61M**

Lot # L328-6

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the GST protein

Cross Reactivity

- All proteins tagged with GST

Host/Isotype/Clone#

Mouse, IgG2a, G2R

Immunogen

Purified GST Protein

Formulation

10mM PBS pH 7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

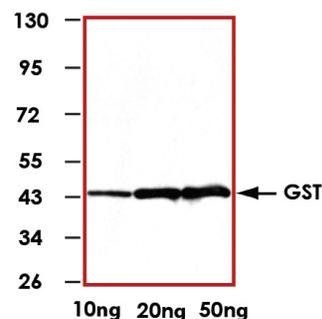
Scientific Background

GST or glutathione S-transferase family of enzymes comprises a long list of cytosolic, mitochondrial, and microsomal proteins that are used to create the so-called 'GST gene fusion system' (1). These proteins are capable of multiple reactions with a multitude of substrates. GST tag has the size of 220 amino acids, which is relatively big compared to other tags like the Myc- or the DYKDDDDK-tag that are quite small. GST can be used to purify and detect proteins of interest.

References

1. Beckett, G.J. and Hayes, J.D. Glutathione S-transferase measurements and liver disease in man. *J. Clin. Biochem. Nutr.* 1987. 2, 1-24.

Sample Data



Representative western blot with Anti-GST (1:1000) using 10 ng, 20 ng, and 50 ng of GST tagged fusion protein.

Anti-GST

Mouse Monoclonal Antibody

Catalog Number

G52-61M

Specific Lot Number

L328-6

Purification

Affinity purified using Protein G

Concentration

1.0 µg/µL

Stability

1yr At -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.