

GST antibody

Cat. No. C15200207

Lot:	002
Size:	500 µg
Type:	Monoclonal
Isotype:	-
Source:	Mouse
Concentration:	1.47 µg/µl

Specificity:	-
Purity:	Protein A purified.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Applications

Applications	Suggested dilution	References
ELISA	1:100,000 – 1:900,000	
Western blotting	1:1,000 – 1:10,000	Fig 1, 2

Target description

GST (Glutathione-S-Transferase) is a protein expression tag commonly used in molecular biology. Anti-GST will react with synthetic construct present in most known GST containing cloning or expression vectors. GST is responsible for the conjugation of reduced glutathione to a wide number of exogenous and endogenous hydrophobic electrophiles. The amino acid sequence GST is highly conserved in most organisms including mammals. GST exists as a 26 kDa homodimer.

Diagenode SA. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders.diagenode@hologic.com
support.diagenode@hologic.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: August, 2024

Results

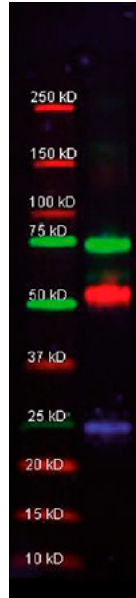


Figure 1: GST antibody western blot results (1/2)

Western Blot of GST antibody. Lane 1: molecular weight marker. Lane 2: Mouseanti- GST monoclonal antibody (blue), Rabbit anti-Transferrin, and Goat-anti-Alpha- 1-Anti-Trypsin were used in a multiplex system to detect target proteins under reducing conditions in albumin depleted human serum with 320 ng of added GST. Load: 1 μ g per lane. Primary antibody: Each primary antibody at 1:1,000 overnight at 4°C. Secondary antibody: DyLight549 Donkey anti-Rabbit IgG (green) DyLight 488 Donkey anti-Mouse IgG (blue), and DyLight 649 Donkey anti-Goat IgG (red) secondary antibody at 1:10,000 for 30 min at RT. Block: MB-070 overnight at 4°C.



Figure 2: GST antibody western blot results (2/2)

Western Blot of GST antibody. Lane 1: GST recombinant protein. Lane 2: lysate of HeLa cells expressing recombinant GST protein. Load: 0.1 μ g per lane. Primary antibody used at 1.0 μ g/ml for 1 h at room temperature. Secondary antibody: IRDye800™ goat anti-mouse secondary antibody at 1:5,000 for 45 min at RT.