

GST monoclonal antibody - Classic

Cat. No. C15200207

Type: Monoclonal

Source: Mouse

Lot #: 001

Size: 500 µg

Concentration: 1.0 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

	Suggested dilution	Results
ELISA	1:100,000 - 1:900,000	
Western blot	1:1,000 - 1:10,000	Figure 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.

Results



Figure 1. GST antibody western blot results

Western Blot of GST antibody. Lane 1: molecular weight marker. Lane 2: Mouse-anti-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 µg 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

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

Diagenode sa. BELGIUM | EUROPE

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

Diagenode Inc. 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: September 10, 2014