

GST polyclonal antibody - Classic

Cat. No. C15430001
Type: Polyclonal
Source: Goat
Lot #: 001
Size: 100 µg
Concentration: 1.0 µg/µl

Specificity: NA

Purity: IgG fraction antibody purified from monospecific antiserum by a multi-step process which includes delipidation, salt fractionation and ion exchange chromatography followed by extensive dialysis

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	1:200 - 1:1,000	
Western blot	1:1000 - 1:5,000	Figure 1

Results

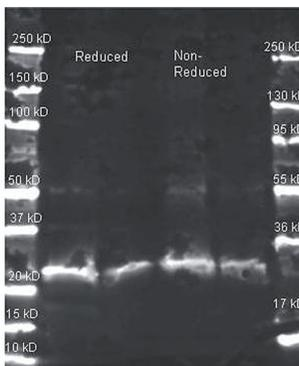


Figure 1. GST antibody western blot results

The Diagenode GST antibody was used to detect GST under reducing and non-reducing conditions. Reduced samples of purified GST contained 4% BME and were boiled for 5 minutes. Samples of ~1 and 0.25 µg of protein per lane were run by SDS-PAGE. Protein was transferred to nitrocellulose and probed with the GST antibody.