
GST Antibody

**Catalog Number
620-0204**

BIO-RAD

Specifications

| | |
|------------------------------|---|
| Source | Affinity-purified mouse IgG _{2a} |
| Formulation | Supplied at 1 mg/ml in phosphate buffered saline, pH 7.4 with 0.05% NaN ₃ . |
| Specificity | The GST-tagged monoclonal antibody recognizes overexpressed GST fusion proteins. |
| Storage and stability | The antibody is shipped on blue ice and should be stored at -20°C for long-term storage. The antibody is stable for 1 year when stored at -20°C (see label for exact expiration). To avoid freeze-thaw cycles, aliquot the antibody. Working solutions can be stored at 4°C for up to 3 months. |
| Volume | 100 µl |

Application

For colorimetric western blotting applications, use at 1:1,000-1:5,000 dilution. For chemiluminescent western blotting applications, use at 1:5,000 dilution.

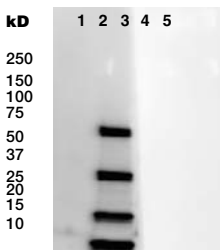


Fig. 1. Chemiluminescent detection of a 26 kD GST protein using the anti-GST antibody at a 1:5,000 dilution. Lane 1: Precision Plus Protein™ standards detected with StrepTactin conjugate; Lanes 2–5: 1,000 ng, 250 ng, 60 ng, and 15 ng of the 26 kD target protein.

StrepTactin is a trademark of Institut für Bioanalytik GmbH.

Expression and purification of GST fusion proteins may require a license under US patent 5,654,176 (assignee: Chemicon International).