## **Profinia**<sup>™</sup> **GST Antibody**

Catalog Number 620-0204



## **Specifications**

**Source** Affinity-purified mouse IgG2a

Formulation Supplied at 1 mg/ml in

phosphate buffered saline, pH 7.4 with 0.05% NaN<sub>2</sub>.

Specificity The GST-tag 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.

## 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.

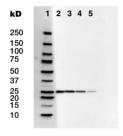


Fig. 1. Chemiluminescent detection of a 26 kD GST protein using the anti-GST antibody at a 1:5,000 dilution. Lane 1: Precision Plus Protein™ standards detected with StrepTactin conjugate, Lanes 2–5: 1,000 ng, 250 ng, 60 ng, and 15 ng of the 26 kD target protein.

