



## **Goat anti-GST (Glutathione-S-Transferase) IgG conjugated to Alkaline Phosphatase**

Product Number    AP-1310  
Antibody            Affinity purified polyclonal Goat anti-GST IgG.  
Amount             100 µg  
Store at             4°C

### Form/Shipping

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Conjugates are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

### Suggested Dilution Range

1:5000-1:50,000 for ELISA and Western blotting

### Handling

Prior to use, reconstitute to 1 mL with ddH<sub>2</sub>O providing a concentration of 100 µg/mL, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

### Buffer

Upon rehydration with 1 mL of ddH<sub>2</sub>O, the product is in 25mM Tris (pH 7.4), 150mM sodium chloride, 5 mM MgCl<sub>2</sub>, 0.12 mM ZnCl<sub>2</sub>, and 2mM sodium azide as a preservative.

### Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months. For extended storage, add an equal volume of glycerol (ACS grade) for a final concentration of 50%, and store at -20 degrees Celsius. Prepare a working dilution daily.

### Country of Origin

USA

### Note

For research use only, not for diagnostic or therapeutic use



**COLUMBIA BIOSCIENCES**