

GST Epitope Tag Antibody

Catalog No. 13-0022
Lot No. 16244001
Pack Size 100 μ l

Type Polyclonal **Host** Rabbit
Mol. Wgt. N/A **Reactivity** N/A
Format Serum **Appl.** IF, WB

Product Description:

Glutathione S-transferase (GST) is a 26 kDa protein from *Schistosoma japonica* and is used as an epitope tag for improving the expression and solubility of GST fusion proteins expressed in *E. coli*. This antibody is useful for detection of GST-tagged proteins on the EpiTitan Histone Peptide array and for detection of GST-tagged fusion proteins by Western blot, including proteins bound to specific peptides in pull-down experiments.

Immunogen:

Purified recombinant GST from *Schistosoma japonica*.

Formulation:

Rabbit serum with 30% glycerol and 0.035% sodium azide.

Storage and Stability:

Stable for 2 years at -20°C from date of receipt.

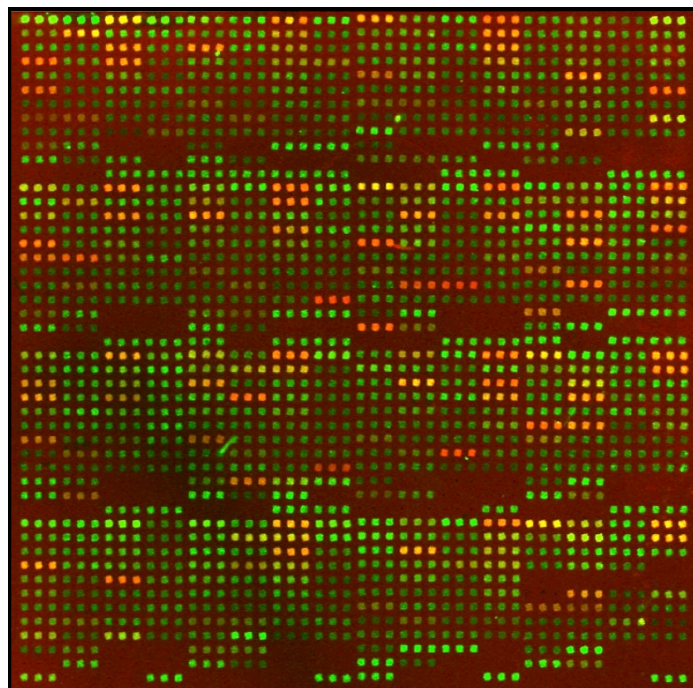
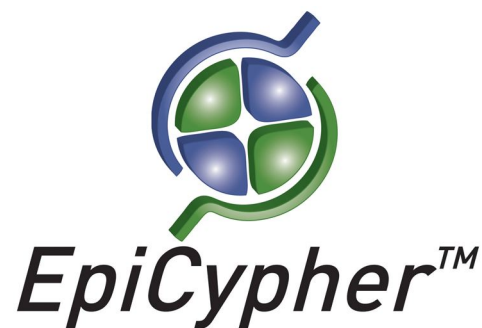
Application Notes:

Recommended Dilution:

IF: 1:750 - 1:2,000
WB: 1:1,000 - 1:5,000

References Using this Product:

Hacker KE et al (2016). *J Biol Chem* DOI: 10.1074/jbc.M116.739375



Immunofluorescence Data: GST Epitope Tag Antibody (1:1,000 dilution) used to label a GST-tagged fusion protein bound to specific peptides on the EpiTitan Histone Peptide Array (Catalog No. 11-2001). The GST antibody was then detected using an AlexaFluor 647-conjugated secondary antibody (red).

Applications Key: ChIP: Chromatin IP; ChIP-seq: Chromatin IP sequencing; E: ELISA; FACS: Flow cytometry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; WB: Western Blotting

Reactivity Key: B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S. cerevisiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish