
GST Tag Antibody [3G10/1B3]

[Catalog No. A26007]

Description

Mouse monoclonal antibody to GST, clone 3G10/1B3

Antibody Type

Monoclonal Antibody

Purification

Purified from hybridoma cell culture supernatant by Protein G chromatography to at least 95% homogeneity as determined by SDS-PAGE.

Immunogen

Recombinant full length protein corresponding to amino acids 1-425 of GST (from *Schistosoma japonicum*, expressed in *E. coli*).

Isotype

IgG1

Size

50 µg

Formulation

Immunogen affinity purified

Specificity

Recognizes Glutathione S-Transferase (GST), a 26-kDa protein encoded by the parasitic helminth *Schistosoma japonicum*.

Storage

Aliquot and store at -20°C for up to 6 months or -80°C for up to 1 year. Avoid repeated freeze/thaw cycles.

Handling Recommendations

For maximum recovery of the products centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

WB: Use at concentration between 0.5 to 2 µg/ml. Detects a band of approximately 30 kDa (predicted molecular weight: 26 kDa); IP: Use at a dilution according to your assay.

Ordering Information:

Products	Size	Cat. No.
GST Tag Antibody [3G10/1B3]	50 µg	A26007-1

This product is for research purposes only. Not intended for use in diagnostic procedures.