

# PRODUCT INFORMATION



## GST-tag Monoclonal Antibody (Clone 2G10B6)

Item No. 37071

### Overview and Properties

<b>Contents:</b>	This vial contains 50, 100, 500 $\mu$ l, or 1 ml of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	Glutathione S-Transferase
<b>Immunogen:</b>	Recombinant <i>S. japonicum</i> GST expressed in <i>E. coli</i>
<b>Species Reactivity:</b>	(+) <i>Schistosoma japonicum</i>
<b>Form:</b>	Liquid
<b>Storage:</b>	-80°C
<b>Stability:</b>	$\geq$ 1 year
<b>Storage Buffer:</b>	0.2 $\mu$ m filtered solution in PBS, with 5% trehalose
<b>Clone:</b>	2G10B6
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG2a
<b>Applications:</b>	ELISA; the recommended starting dilution is 1:1,000-1:2,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Description

GST-tag Monoclonal Antibody (Clone 2G10B6) is a probe for the immunochemical detection of GST tags on recombinant proteins. Recombinant proteins are commonly labeled with affinity tags, such as GST, to facilitate both their detection and purification.<sup>1</sup> GST tags are commonly utilized to increase protein solubility and enhance expression efficiency. Cayman's GST-tag Monoclonal Antibody (Clone 2G10B6) can be used for ELISA. It recognizes proteins containing GST tags, proteins containing GST- and His-tags, and GST with an HA-tag at either the N- or C-terminus.

### Reference

1. Zhao, X., Li, G., and Liang, S. Several affinity tags commonly used in chromatographic purification. *J. Anal. Methods Chem.* **581093** (2013).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/05/2022

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM