

Product Information



Glutathione S-Transferase Polyclonal FITC Antibody

Item No. 10197

Contents:	This vial contains 100 µg of fluorescein-labeled IgG in 500 µl PBS, pH 7.2, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Synonym:	GST
Antigen:	Purified glutathione S-transferase (<i>S. japonicum</i>)
Host:	Rabbit
Cross Reactivity:	(+) GST and GST fusion proteins; other species not tested
Stability:	≥1 year at 4°C
Applications:	The recommended starting dilution for western blot and immunofluorescence is 1:200-1:400 (1.0-0.5 µg/ml). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Cayman's Glutathione S-Transferase (GST) Polyclonal FITC Antibody may be used to monitor expression of recombinant proteins tagged with GST (GST fusion proteins). A recombinant GST-tagged protein is expected to migrate 26 kDa slower than the non-tagged version in reducing conditions on SDS-PAGE. In general, GSTs are ubiquitous multifunctional enzymes, which play a key role in cellular detoxification. Researchers routinely utilize the enzyme as a method to simplify the purification of recombinant proteins.¹

Reference

1. Boyland, E. and Chasseaud, L.F. The role of glutathione and glutathione S-transferases in mercapturic acid biosynthesis. *Adv. Enzymol.* **32**, 173-219 (1969).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10197

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/17/2011