

PRODUCT INFORMATION



GST-tag Polyclonal Antibody (FITC)

Item No. 10197

Overview and Properties

Contents:	This vial contains 500 µl of fluorescein-labeled polyclonal antibody.
Synonyms:	Glutathione S-Transferase
Immunogen:	Purified glutathione S-transferase (<i>S. japonicum</i>)
Cross Reactivity:	(+) GST and GST fusion proteins
Species Reactivity:	(+) <i>S. japonicum</i> ; other species not tested
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunofluorescence (IF) and Western blot (WB); the recommended starting dilution for IF and WB is 1:200-1:400 (1.0-0.5 µg/ml). Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

GST-tag Polyclonal Antibody (FITC) 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.¹ GST tags are commonly utilized to increase protein solubility and enhance expression efficiency. Cayman's GST-tag Polyclonal Antibody (FITC) is composed of a GST-tag polyclonal antibody conjugated to fluorescein isothiocyanate (FITC) and can be used for immunofluorescence (IF) and Western blot (WB) applications. It recognizes GST and proteins containing GST tags.

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, 07/06/2023

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