

Glutathione S-transferase monoclonal antibody (8-326) (biotin conjugate)

Epitope tags provide a method to localize gene products in a variety of cell types, study the topology of proteins and protein complexes, identify associated proteins, and characterize newly identified, low abundance or poorly immunogenic proteins when protein specific antibodies are not available. The GST (Glutathione-S-Transferase) is a 26kDa protein encoded by the parasite *Schistosoma japonicum*, easily purified by affinity chromatography on immobilized glutathione. GST is utilized as a fusion protein with foreign proteins in a range of prokaryotic expression vectors, including the pGEX family of vectors. Antibodies against GST will detect the expressed GST protein or a GST-fusion protein, and can be utilized for immuno-purification of GST-fusion proteins.

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GST
Application	IP, WB
Clone	8-326
Formulation	Liquid. In PBS containing a proprietary stabilizer and 0.02% sodium azide.
Host	Mouse
Immunogen	Full-length recombinant human Glutathione S-transferase.
Isotype	IgG
Purity Detail	Purified
Species Reactivity	Species independent
UniProt ID	P08263
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com