

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](#).

Ordering Information

[Order Online »](#)

ENZ-ABS280-0050	50µl
-----------------	------

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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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