

GST-P polyclonal antibody

The GST superfamily of enzymes play an important role in detoxification of xenobiotics by catalyzing the conjugation of many hydrophobic and electrophilic compounds with reduced glutathione. There are four main GST classes; alpha, mu, pi, and tau. GST-P is a member of the pi class of GSTs, and its expression is specifically increased in the early stages of rat hepatocarcinogenesis. Due to its high affinity for glutathione, GST is frequently utilized as an approximately 26 kDa fusion tag for the expression and purification of recombinant proteins.

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

Citations: 5

[View Online »](#)

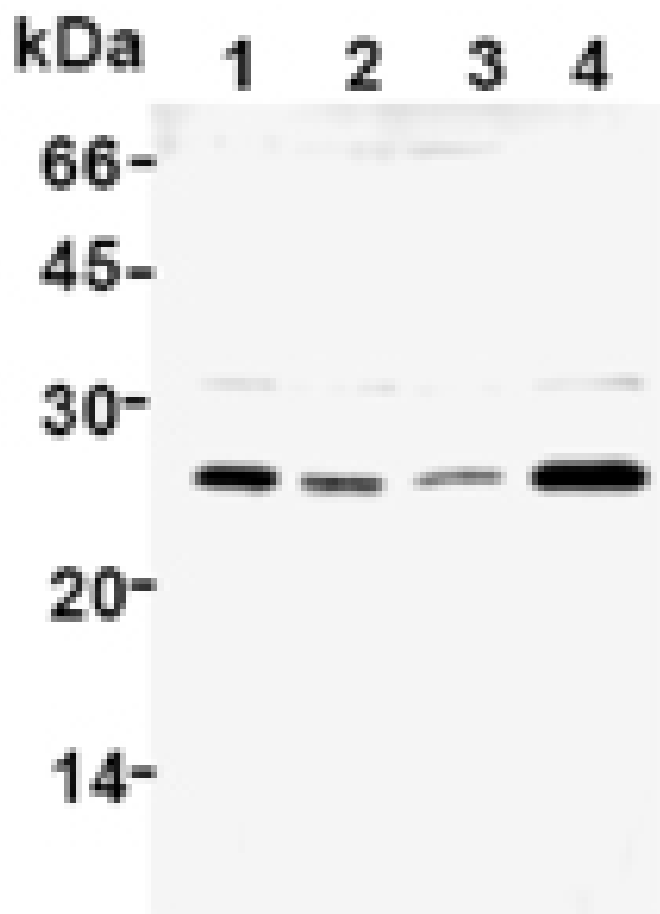
Ordering Information

[Order Online »](#)

ADI-MSA-101-E	100µl
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of mouse kidney (1), WR19L (2), PC12 (3) and Jurkat (4) using GST-P pAb.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name GST-π, Glutathione S-transferase P, GSTP1

Application IHC, WB

Application Notes Detects a band of ~27kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol.

Host Rabbit

Immunogen Native rat liver GST-P.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:1,000-1:2,000)Western Blot (1:1,000-1:2,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P04906

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](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-
be@enzolifesciences.com](mailto:info-be@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)