

# GST- $\pi$ 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: 24

[View Online »](#)

## Ordering Information

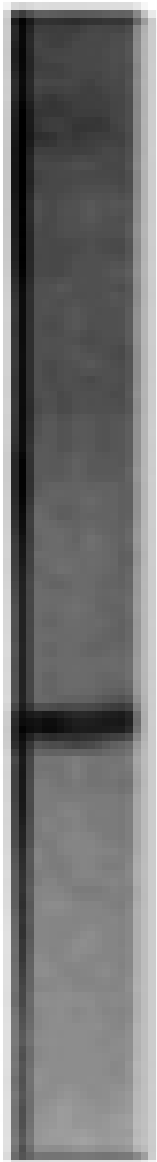
[Order Online »](#)

ADI-MSA-102-E	100 $\mu$ l
---------------	-------------

## Manuals, SDS & CofA

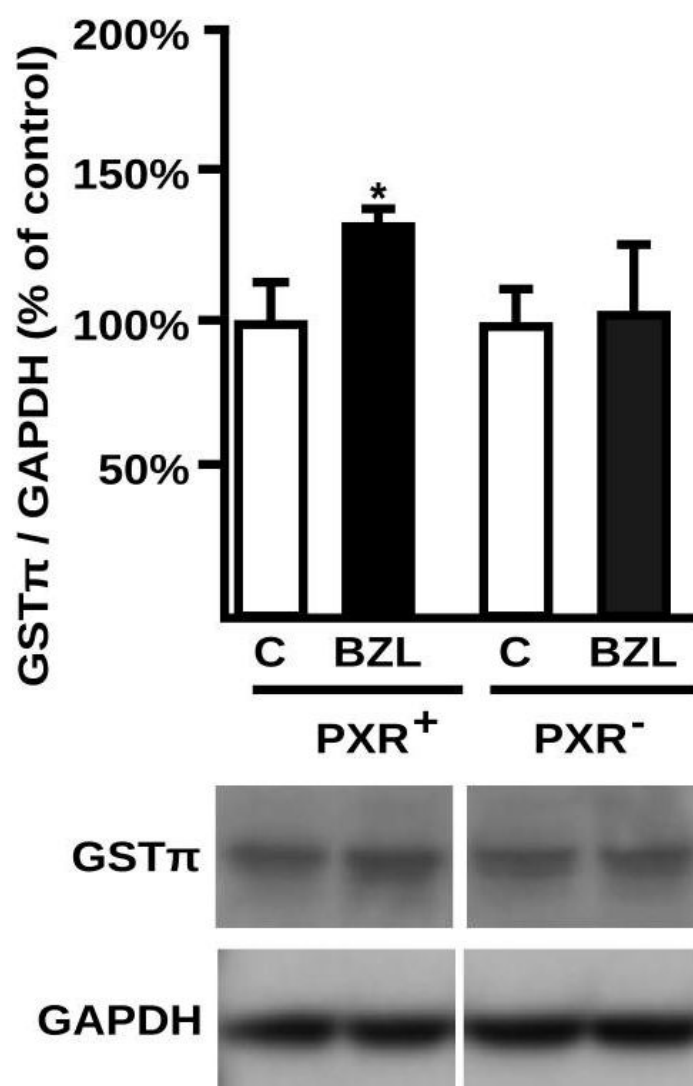
[View Online »](#)





Western blot analysis of mouse liver tissue extract, probed with GST-p pAb.

**D**



Effect of PXR knock down on BZL mediated P-gp, MRP2, CYP3A4 and GSTπ induction. P-gp (panel A), MRP2 (panel B), CYP3A4 (panel C) and GSTπ (panel D) levels were estimated by western blotting in lysates from HepG2 cells transfected either with 100 nM Control siRNA-A (PXR<sup>+</sup>) or 100 nM PXR siRNA (h) (PXR<sup>-</sup>) and exposed to BZL (200 μM, 48 h) or vehicle (C). Equal amounts of total protein (7 μg) were loaded in the gels. O.D. from each protein was related to GAPDH O.D. Uniformity of loading and transfer from gel to PVDF membrane was also controlled with Ponceau S. Typical western blot detections from each group are shown at the bottom of bar graphics. The results (% of each control) are expressed as mean ± S.D. (n = 3). \*Significantly different from C, p<0.05.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Regulation of biotransformation systems and ABC transporters by benznidazole in HepG2 cells: involvement of pregnane X-receptor. *PLoS Negl Trop Dis* (2012)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**Application** IHC, WB

**Application Notes** Detects a band of ~26kDa by Western blot.

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol.

**Host** Rabbit

**Immunogen** Human placental GST- $\pi$ .

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:500-1:1,000)Western Blot (1:500-1:1,000)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** P09211

**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-  
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)