



**EAGLE**  
BIOSCIENCES

**Anti-Rat GST Ya Antiserum**

Catalog Number: GST00-A100

Aliquot: 100  $\mu$ L

Storage: -70°C

**PRODUCT SPECIFICATION SHEET**

<b>DESCRIPTION:</b>	Goat antisera specific for Rat Glutathione S Transferase Ya. This antibody does not cross react with the mu or pi class of GST.
<b>CONCENTRATION:</b>	Not Determined
<b>STORAGE:</b>	-70°C. <b>AVOID REPEATED FREEZE THAWS.</b>
<b>STORAGE BUFFER:</b>	None
<b>IMMUNOGEN:</b>	Affinity Purified Rat Liver GST Ya
<b>PURIFICATION:</b>	None
<b>APPLICATIONS:</b>	Western Blotting: 1:1000 This antiserum yields a positive result with 0.1 $\mu$ g of Purified Rat GST Ya when using a 1:1000 dilution of GS 09 and a 1:5000 dilution of alkaline phosphatase-conjugated anti-goat secondary antibody.
<b>SPECIES SPECIFICITY:</b>	Rat and Human.
<b>REFERENCE:</b>	Holmes, W., <i>et. al.</i> , <i>Br. J. Cancer</i> , <b>62</b> :209-212 (1990)

*Product Developed and Manufactured in the USA*

---

EAGLE BIOSCIENCES, INC.  
20A NW Blvd., Suite 112  
Nashua, NH 03063

Phone: 617-419-2019  
Fax: 617-419-1110  
Email: [info@EagleBio.com](mailto:info@EagleBio.com)  
[www.EagleBio.com](http://www.EagleBio.com)