



EAGLE
BIOSCIENCES

Anti-Phosphotyrosine IgG

Catalog Number: PHT00-A250

Aliquot: 250 μ g

Storage: -70°C

PRODUCT SPECIFICATION SHEET

DESCRIPTION:	This goat polyclonal antibody is specific for phosphotyrosine
MOLECULAR WEIGHT:	The MW of bands on a Western blot is dependent upon the protein being probed.
CONCENTRATION:	0.5 mg/mL as determined using BCA assay with BSA as a reference standards.
STORAGE BUFFER:	PBS
STORAGE:	-70°C. AVOID REPEATED FREEZE THAWS.
IMMUNOGEN:	Phosphotyrosine
CLASS:	IgG
PURIFICATION:	Ammonium Sulfate Fractionation
APPLICATIONS:	Suitable for use in Western blotting, ELISA and immuno-precipitation applications.
SPECIES SPECIFICITY:	Not species specific. May not exhibit utility when using in a goat experimental model.

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
www.EagleBio.com