



018PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

AP Conjugated Rabbit α -Human IgG

Affinity Purified Secondary Antibody

(Cat. # 786-R49)



think proteins! think G-Biosciences www.GBiosciences.com

DESCRIPTION

Affinity purified rabbit α -human IgG antibody-Alkaline phosphatase conjugate.

ITEM(S) SUPPLIED

Cat. #	Description	Concentration	Final Volume
786-R49	AP labeled rabbit α -human IgG	1mg/ml	1ml

STORAGE

The conjugate should be stored at 4°C and is stable for up to 1 year. DO NOT FREEZE.

PURIFICATION:

The antibody is isolated from antisera by immunoaffinity chromatography using human IgG coupled to agarose beads followed by solid phase adsorption(s) to remove unwanted reactivities.

FORMAT

A sterile filtered liquid. The storage buffer is 50mM Tris.HCl, 150mM NaCl, 1mM MgCl₂, 0.1mM ZnCl₂, 50% glycerol (v/v) at pH 8.0. This is supplemented with 10mg/ml Bovine serum albumin (BSA) as a stabilizer and 0.1% (w/v) sodium azide as a preservative.

APPLICATIONS/RECOMMENDED WORKING DILUTIONS

Application	786-R49
ELISA	1:2,000-1:10,000
Immunohistochemistry	1:200-1:1,000
Western blot	1:500-1:2,500

NOTE: Optimal working dilution of AP-conjugate for the specific application must be determined by the end user to obtain the best conditions. Working diluted solution should be prepared immediately before use and unused diluted solution should be discarded.

RELATED PRODUCTS

Download our Western Blotting and Assay Development Handbooks.



<http://info.gbiosciences.com/complete-western-blot-handbook--selection-guide>

<http://info.gbiosciences.com/complete-assay-development-handbook/>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 10/3/2012 CMH



www.GBiosciences.com