



017PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# AP Conjugated Affinity Purified Antibody

## Secondary Antibody

(Cat. # 786-R43 to 786-R47, 786-R49)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## DESCRIPTION

Affinity purified antibody-Alkaline phosphatase conjugate (as labeled).

## ITEM(S) SUPPLIED

Cat. #	Description	Concentration	Final Volume
786-R43	AP labeled goat $\alpha$ -mouse IgG	0.6mg/ml	1ml
786-R44	AP labeled goat $\alpha$ -rabbit IgG	1mg/ml	1ml
786-R45	AP labeled goat $\alpha$ -rat IgG	0.6mg/ml	1ml
786-R46	AP labeled goat $\alpha$ -human IgG	0.6mg/ml	1ml
786-R47	AP labeled rabbit $\alpha$ -goat IgG	0.6mg/ml	1ml
786-R49	AP labeled rabbit $\alpha$ -human IgG	<a href="#">Download separate protocol</a>	

## STORAGE

The conjugate should be stored at 4°C and is stable for up to 1 year. After reconstitution the product is stable for several weeks at 2-8°C as an undiluted liquid. For extended storage after reconstitution, add an equal volume of glycerol to make a final concentration of 50% glycerol (ACS grade or better) followed by storage at -20°C.

**NOTE:** *The concentration of antibody (and buffers) will decrease to one-half of the original after the addition of glycerol. Avoid repeated freeze/thaw cycles*

## PURIFICATION:

The antibody is isolated from antisera by immunoaffinity chromatography using antigen coupled to sepharose beads.

## FORMAT

Lyophilized in a buffer containing 0.01M Sodium Phosphate, 0.25M NaCl, pH 7.6 with 15mg/ml BSA and 0.01% Thimerosal.

## RECONSTITUTION

Reconstitute with sterile distilled water to match the Final Volume indicated in the table above.

## APPLICATIONS/RECOMMENDED WORKING DILUTIONS:

- EIA and Western blots: 1:5,000-1:100,000.
- Immunohistochemistry: 1:500-1:5,000.

**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.*

## WARNING

**NOTE:** Use of sodium azide, an inhibitor of phosphatases will substantially inhibit the enzyme activity of AP.

## 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](http://www.GBiosciences.com) or contact us.

Last saved: 10/3/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)