

BIO-RAD

Affinity Purified
Goat anti-human IgG (γ)
Catalog No. 172-1052
Control No. 26274

Contents: 1 mg Antibody

Reconstitution: For most purposes, the lyophilized contents of the vial should be reconstituted with 1 ml of sterile, 40% glycerol. However, if the antibody is to be labeled with an enzyme or other compound, it is recommended that the antibody be rehydrated and annealed by the following procedure.

Add 0.1 ml of **0.01M** acetic acid to the vial and rotate the vial until the lyophile is dissolved.

Add 0.1 ml of freshly prepared **0.2M NaCl** in **0.02M Na₂CO₃**. Mix rapidly to obtain a clear or opalescent solution. The **pH** should be approximately 9.5.

Replace the stopper and place the vial in a 400 ml water bath preheated to **55-60°C**. Place the bath on the bench and allow it to slowly cool to room temperature.

Add 0.8 ml of phosphate buffered saline or other desired buffer. The final **pH** should be approximately 7.5.

Storage: The unopened vial of lyophilized antibody can be stored at 4°C. Antibody reconstituted with glycerol, as in the first procedure above, can also be stored at 4°C. Rehydrated antibody should be stored frozen, but repeated freeze-thaw cycles should be avoided.

Source of Antibodies: These antibodies were isolated from a pool of sera from goats immunized with purified **IgG** from pooled normal human sera.

This product is intended for research use only. It is not intended for clinical diagnostic purposes.

Bio-Rad Laboratories, 32nd & Griffin Ave., Richmond, CA 94804