

BIO-RAD

**Affinity Purified
Goat anti-Human IgE (ϵ)
Catalog No. 172-1061**

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 a purified IgE kappa myeloma protein and the Fc portion of an IgE lambda myeloma protein.

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

86-0282 586

Bio-Rad Laboratories, 32nd & Griffin Ave., Richmond, CA 94904