

BIO-RAD

***EIA Grade Affinity Purified Antibody
Goat Anti-Rabbit IgG (H + L)
Human IgG Adsorbed***

Catalog No. 172-1053

- Contents:** 2 mg Antibody
- Reconstitution:** For most purposes, the lyophilized contents of the vial should be reconstituted with 2 ml of sterile, 40% glycerol.
- Storage:** The unopened ~~vial of lyophilized~~ antibody can be stored at **4°C**. Antibody reconstituted with glycerol can be stored at 4°C or -20%.
- source of Antibodies:** These antibodies were isolated from a pool of sera from goats immunized with purified IgG from pooled normal rabbit sera

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