
Immun-Star™
Goat Anti-Rabbit
(GAR)-HRP Conjugate

Catalog #
170-5046

BIO-RAD

Contents — 2 ml lyophilized conjugate

Antibody Concentration — 0.8 mg/ml
after reconstitution

Buffer composition after reconstitution
— 0.01 M sodium phosphate, 0.25 M NaCl,
pH 7.6, 0.01% Thimerosal, 15 mg/ml BSA.

Antibody preparation — The antibody has
been isolated from antisera by immunoaffinity
chromatography using antigen coupled to
agarose beads.

Antibody specificity — This antibody
reacts with the heavy chains of rabbit IgG
and with the light chains common to most
rabbit immunoglobulins. No antibody was
detected against non-immunoglobulin serum
proteins; however, this antibody may cross-
react with other species.

Shelf life — One year at 4°C

Reconstitution — Restore with 2 ml distilled water. Centrifuge product if not completely clear after standing for 1–2 hours at room temperature.

Recommended working dilution — 1:20,000–1:100,000 for Immun-Star HRP chemiluminescent kit.

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Drive, Hercules, CA 94547 USA

510-741-1000

10000090 Rev A