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

Catalog Number
170-6401

Specifications

Contents	1.0 ml antibody solution.
Preparation of antibody	Goat anti-rabbit IgG (H+L) is prepared from antisera raised in goats immunized with purified rabbit IgG. Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies that cross-react with human immunoglobulin.
Biotinylation of antibody	Antibody is biotinylated using a modification of the procedure described by Guesdon et al. ¹
Buffer	10 mM Tris, 150 mM sodium chloride, pH 7.5. Contains 1.0% bovine serum albumin as a stabilizer and 0.1% sodium azide as a bacteriostat.

Storage conditions

This reagent is shipped frozen on dry ice and can be stored at -20 °C prior to opening. Once thawed, store at 4 °C. Repeated freeze-thaw cycles should be avoided.

Shelf life

1 year at 4 °C.

Recommended working dilution

1:3,000 as determined by western blot.

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

Reference

1. Guesdon, J. L., Ternyck, T. and Avrameas, S., *J. of Histochemistry and Cytochemistry*, **27**:8, 1131–1139 (1979).