
Blotting Grade Affinity Purified Goat Anti-Mouse IgG (H+L) Alkaline Phosphatase Conjugate

Catalog Number
170-6520

Bio-Rad Laboratories, 2000 Alfred Nobel Dr., Hercules, CA 94547

LIT420 Rev D

BIO-RAD

Specifications

Contents 1 ml

Buffer 10 mM Tris, 150 mM sodium chloride, 1.0 mM magnesium chloride, pH 8.0. Contains 1.0% bovine serum albumin and 0.1% sodium azide.

Preparation of antibody Goat Anti-Mouse IgG (H+L) is prepared from antisera raised in goats immunized with purified mouse IgG. Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies which cross react with human immunoglobulin.

Preparation of conjugate Antibody is conjugated to alkaline phosphate using a conjugate modification of the

one step glutaraldehyde procedure described by Engvall and Perlmann.¹

Storage Goat Anti-Mouse IgG (H+L) AP conjugate will arrive frozen. It may be stored frozen until needed. Once thawed, store at 4 °C. Repeated freeze-thaw cycles should be avoided.

Shelf life One year, at 4 °C.

Recommended working dilution 1:3,000

Reference

1. Engvall, E. and Perlmann, P., *Immunochemistry*, 8, 871 (1971).