

Anti-Human IgG (Goat), Biotin-Labeled	NEF803001EA
Anti-Rabbit IgG (Goat), Biotin-Labeled	NEF813001EA
Anti-Mouse IgG (Goat), Biotin-Labeled	NEF823001EA

STORAGE CONDITIONS:

These products are supplied lyophilized and should be stored at 4 C until rehydrated. Bovine serum albumin is added as a protein stabilizer.

PRODUCT DESCRIPTION:

These antibodies were affinity purified from goat serum immunized with purified human, rabbit or mouse IgG, and labeled with a biotin analog having a seven atom linker consisting of 6-aminocaproic acid and the 6 atom side chain of biotin. Each vial contains 0.5 mg of affinity purified antibody. These products are tested for specificity by immunoelectrophoresis, gel diffusion and ELISA techniques. No antibody is detectable to non-immunoglobulin serum components.

REHYDRATION AND STORAGE:

Two methods for rehydration and storage are recommended to meet most needs:

Procedure A: uses 50% glycerol and allows 4 C or lower storage for several months.

Procedure B: employs water for procedures in which glycerol may deter maximum function or where a frozen product is preferred for long term storage.

Procedure A:

Mix 1 ml of glycerol with 1 ml of water in a test tube. Transfer 1 ml of the 50% glycerol solution to the product vial. Rotate the vial until the lyophilized pellet is totally dissolved. This product may be stored either in the refrigerator or in the freezer as desired.

Procedure B:

Rehydrate with 1 ml of reagent quality water. This product may be stored for up to one week refrigerated; thereafter, it should be stored frozen.

PERFORMANCE CHARACTERISTICS:

Biotin-labeled affinity purified antibodies are excellent for both histochemical/cytochemical staining procedures and for use in enzyme amplified immunoassays such as microplate and membrane ELISA. Serial dilutions of all antibody reagents should be performed to determine the best working concentrations under a given set of assay conditions. Suggested starting concentrations are as follows:

Microplate and Membrane Immunoassays: 1:1000
Histo/Cytochemical Procedures 1:50 - 1:500

HAZARD WARNING:

This product is considered to be non-hazardous. Although the product is classified as nonhazardous, we strongly recommend using prudent laboratory practices: Avoid unnecessary contact, use of gloves, eye protection, lab coats, etc. when using this or any laboratory reagent.

Limited Warranty: This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.