

Caution: For Laboratory Use. A product for research purposes only

Alkaline Phosphatase Labeled Affinity Purified Antibody to Rabbit IgG (H+L) [Goat]

Product Number: NEF814001EA

Alkaline Phosphatase Labeled Affinity Purified Antibody to Rabbit IgG (H+L) [Goat]

LOT NO:

SIZE 1.0 mL

DESCRIPTION Affinity purified antibody, isolated from a pool of serum from goats immunized with purified rabbit IgG was labeled with phosphatase using a modified glutaraldehyde procedure.

FORM/STORAGE Liquid. Store refrigerated at 2 – 8°C.

CONTENT Each vial contains 1.0 mL at 1.0 mg/mL of affinity purified antibody from unlabeled antibody

LOT NO

E/P RATIO

The calculated enzyme/antibody protein ratio:

2.40: 1

SPECIFICITY

Tested by immunoelectrophoresis, gel diffusion and ELISA techniques as applicable. This product reacts specifically with rabbit IgG, heavy and light chains. No antibody is detectable to non-immunoglobulin serum components.

SUGGESTED WORKING DILUTIONS

Optimal working concentration should be determined experimentally. Prepare working dilution in TBS or other buffer such as blocking solution. These buffers are not recommended for long term storage. Suggested starting dilutions are as follows. **Note:** In many cases, the antibody may be diluted more than indicated.

Microwell:

1 / 2,000 – 1 / 10,000 0.5 µg/mL – 0.1 µg/mL

Blotting:

1 / 10,000 – 1 / 40,000 0.1 µg/mL – 0.025 µg/mL

Histo / Cytochemical Procedures:

1/400 – 1/1,000 2.5 µg/mL – 1.0 µg/mL

PRODUCT SAFETY AND HANDLING

This product is considered non-hazardous as defined by The Hazard Communication Standard (29 CFR 1910.1200). Avoid contact with skin and eyes. In case of contact or spillage, clean with copious amounts of water. Dispose via sanitary sewer.

STABILIZERS AND PRESERVATIVES

Goat serum and bovine serum albumin added as protein stabilizer. No preservatives added. Non-sterile.

STABILITY

Product is stable for one year from date of receipt when stored properly.

COUNTRY OF ORIGIN OF ANTISERA United States

DATE OF MANUFACTURE:

PerkinElmer, Inc.
549 Albany Street
Boston, MA 02118 USA
P: (800) 762-4000 or (+1) 203-925-4602
www.perkinelmer.com/nenradiochemicals
For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs
Copyright ©2010, PerkinElmer, Inc. All rights reserved.
PerkinElmer® is a registered trademark of PerkinElmer, Inc.
All other trademarks are the property of their respective owners.

