



CERTIFICATE OF ANALYSIS

Product ANTI-RAT IgG (H+L)
 AFFINITY PURIFIED

Catalog AI-4000

Amount 1.5 mg IgG

Concentration 3.0 mg/ml

Lot No. ZC0331

Produced in rabbit

Solution 10 mM phosphate, pH 7.8, 0.15 M NaCl, 0.08% sodium azide. Sucrose has been added for stabilization.

Storage Refrigerate; for long-term storage, aliquot and store frozen.

Antibody Activity (by solid phase IgG binding assay) 96%

Solid-phase enzyme immunoassay of cross-reactivity:

<u>Protein</u>	<u>Percent Cross-reactivity</u>
hamster IgG	12%
bovine IgG, guinea pig IgG, mouse IgM, Rhesus monkey IgG	6%
cat IgG, horse IgG, mouse IgG, swine IgG	3%
goat IgG, sheep IgG, swine IgM	1%
bovine IgM, chicken IgG, dog IgG, dog IgM, goat IgM, guinea pig IgM, horse IgM, human IgG, human IgM, rat IgM, sheep IgM	<1%

If this antibody is to be used in tissues which may contain cross-reacting endogenous immunoglobulins, dilution of this antibody may be made in buffers containing 2 % normal serum from the same species as the tissue.

The recommended concentration range for use is 5-20 µg/ml.



CERTIFICATE OF ANALYSIS