
Non-Immune Mouse IgG Secondary Antibody

(Base Catalog No. A16000)

Description

Non-Immune Mouse IgG has wide applications in the routine immunocytochemical detection of mouse monoclonals and can be used with anti-mouse Ig bridging antibody, raised in any appropriate species, such as rabbit, sheep or goat. It can be used in both light and electron microscopes. Dilutions of 1/500 - 1/5000 have produced excellent results using a wide range of primary monoclonal antibodies. The optimum dilution should be determined for each primary monoclonal antibody and anti-mouse bridging antibody used.

Purification

Purified mouse IgG conjugated to Horseradish Peroxidase (HRP) - liquid

Isotype

IgG

Concentration

1 mg/ml , 0.5 ml

Formulation

Phosphate buffered saline. 0.01% Thiomersal ; HRP Stabiliser

Storage

Store at 4°C ; DO NOT FREEZE; This product should be stored undiluted

Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

ELISA, ChIP

Ordering Information:

Products	Size	Cat. No.
Non-Immune Mouse IgG Secondary Antibody	0.5 ml	A16000-1

This product is for research purposes only. Not intended for use in diagnostic procedures.