
Non-Immune Rabbit IgG Secondary Antibody

(Catalog No. A16001)

Description

Purified Rabbit IgG

Purification

Purified polyclonal rabbit IgG - liquid

Isotype

Polyclonal IgG

Concentration

0.5 ml

Formulation

Phosphate buffered saline. 0.09% Sodium Azide

Specificity

Rabbit

Storage

Store at 4°C or at -20°C if preferred ;Storage in frost-free freezers is not recommended;This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody.

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 Rabbit IgG Secondary Antibody	0.5 ml	A16001-1

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