

Datasheet for Rabbit Anti-Chicken IgY Fab

Catalog number: mAb-02616-100

Background: Secondary antibodies are affinity-purified antibodies which will work with target-specific primary antibody in the detection, sorting or purification of its specified target. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies. Most commonly, secondary antibodies are generated by immunizing the host animal (different from host species of primary antibody) with a pooled population of normal immunoglobulins from the host species of primary antibody and can be further purified and modified (i.e. antibody fragmentation, label conjugation, etc.) to ensure well-characterized specificity to corresponding normal immunoglobulins.

■ Product Information

Symbol: Rabbit Anti-Chicken

Alias:

Cross reactivity:

Expected Protein Weight:

Speices:

Source: Rabbit

Isotype: IgG

Purity: Affinity purification

Storage: Store at -20°C. Avoid freeze / thaw cycles. | Buffer: PBS with 0.05% proclin300, 0.05% BSA, 50% glycerol, pH7.3.

■ Immunogen Information

Gene ID:

Swiss:

Immunogen: Chicken IgY

■ Recommended Dilutions

WB 1:2000 - 1:20000

Citation of product: If use of this item results in a publication, please use this information: Rabbit Anti-Chicken IgY Fab(mAb-02616-100; Genecopoeia, Inc, Rockville, MD).

GeneCopoeia Products are for Research Use Only
9620 Medical Center Drive, #101 Rockville, MD 20850 USA
Tel: 301-762-0888: Fax: 301-762-3888
Email: support@genecopoeia.com; Web: www.genecopoeia.com

Copyright © 2024 GeneCopoeia, Inc.