

Anti-Rabbit IgG Fc Antibody (14H9H10)[HRP], mAb, Mouse

Abbreviated name-1: Anti-Rabbit IgG Fc Monoclonal Secondary Antibody (Min X Hu, Ms, Rt, Sh, Bv, Gt, Camel)

PRODUCT INFORMATION

Description

Generally, goat anti-rabbit or donkey anti-rabbit IgG polyclonal antibodies are used as secondary antibodies to detect rabbit primary antibodies. However, since polyclonal antibodies can recognize many epitopes, the background and the cross reactivity in assays is usually high, making results hard to interpret.

To minimize the background, cross reactivity, and improve the lot-to-lot consistency, a unique **GenScript Anti-Rabbit IgG Fc Antibody (14H9H10)[HRP], mAb, Mouse** has been developed to be used as the secondary antibody. Simply substitute conventional HRP or fluorescent reagents for unparalleled clarity in results and high quality images in immunoprecipitation, Western Blot, ELISA, immunohistochemistry and immunofluorescence.

Specificity

GenScript Anti-Rabbit IgG Fc Antibody (14H9H10)[HRP], mAb, Mouse reacts with the Fc portion of rabbit IgG but not with the Fab portion of rabbit IgG. The monoclonal antibody has no cross-reactivity with human, mouse, rat, sheep, bovine, goat or camel immunoglobulins.

Concentration

0.5 mg/ml, lyophilized with PBS, pH 7.4, containing 0.02% sodium azide. (It could be customized for bulk order)

Storage

GenScript Anti-Rabbit IgG Fc Antibody (14H9H10)[HRP], mAb, Mouse should be stored at -20°C. This product is stable for 1 year upon receipt, when handled and stored as instructed.

Applications

Working concentrations for specific applications should be determined by the investigator. The appropriate concentrations may be affected by primary antibody affinity, antigen concentration, the sensitivity of the method of detection, temperature, the length of the incubations, and other factors. The suitability of this antibody for applications other than those listed below has not been determined. The following concentration ranges are recommended starting points for this product.

Western Blotting: 1:20,000-1:40000

ELISA: 1:20,000-1:40,000

Immunohistochemistry: user-optimized

Optimal working dilutions must be determined by the end user.

Cat. No.: A01856-40

Host: Mouse

Size: 40 µg

Immunogen: Rabbit IgG(H&L)

Ig Subclass: IgG1, κ

Clone: 14H9H10

Purification: Protein A affinity column

Conjugation: HRP

Version: 01/03/2017

** For non-clinical research use only. **

860 Centennial Ave., Piscataway, NJ 08854, USA