

## Anti-Human IgG Fc Antibody (50B4A9)[HRP], mAb, Mouse

**Abbreviated name-1: Anti-Human IgG Fc Monoclonal Secondary Antibody (Min X Ms, Rt, Rb, Ch, Gt)**

### PRODUCT INFORMATION

#### Description

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

To minimize the background and cross reactivity and improve the lot-to-lot consistency, a unique **GenScript Anti-Human IgG Fc Antibody (50B4A9)[HRP], mAb, Mouse** has been developed to be used as a secondary antibody. The superiority of this antibody is that it can be used as an idiotype antibody to detect drug metabolism *in vivo*.

#### Specificity

**GenScript Anti-Human IgG Fc Antibody (50B4A9)[HRP], mAb, Mouse** reacts with the Fc portion of human IgG but not with the Fab portion of human IgG. This monoclonal antibody has no cross-reactivity with mouse, rat, rabbit, chicken or goat 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-Human IgG Fc Antibody (50B4A9)[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: 1:5,000-1:20,000

ELISA: 1:5,000-1:20,000

Immunohistochemistry: user-optimized

Optimal working dilutions must be determined by the end user.

**Cat. No.:** A01854-40

**Host:** Mouse

**Size:** 40 µg

**Immunogen:** Human IgG(H&L)

**Ig Subclass:** IgG2b,k

**Clone:** 50B4A9

**Purification:** Protein A affinity column

**Conjugation:** HRP

**Version:** 01/03/2017

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

**860 Centennial Ave., Piscataway, NJ 08854, USA**