

## Anti-Human IgG Fab Antibody (12H3C4A6) [HRP], mAb, Mouse

**Abbreviated name-1: Anti-Human IgG Fab 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 antibodies to detect human IgG (H+L). However, since polyclonal antibodies can recognize many epitopes, the background and the cross reactivity in assays is usually high, making results hard to interpret. As a mouse monoclonal antibody, the unique **GenScript Anti-Human IgG Fab Antibody (12H3C4A6)[HRP], mAb, Mouse** has a high specificity to the human IgG Fab portion and can be used to detect human antibodies which include the human IgG Fab portion.

#### Specificity

**GenScript Anti-Human IgG Fab Antibody (12H3C4A6)[HRP], mAb, Mouse** reacts with the Fab portion of human IgG but not with the Fc portion of human IgG. The 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 Fab Antibody (12H3C4A6)[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.:** A01855-40

**Host:** Mouse

**Size:** 40 µg

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

**Ig Subclass:** IgG1, $\kappa$

**Clone:** 12H3C4A6

**Purification:** Protein A affinity column

**Conjugation:** HRP

**Version:** 01/03/2017

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

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