

**TRITC Anti-human CD40 Antibody \*G28.5\***Catalog number: 104011J0, 104011J1  
Unit size: 100 tests, 500 tests**Product Details**

|                    |                                                                             |
|--------------------|-----------------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving                                                    |
| Concentration      | 0.1 mg/mL                                                                   |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

|                    |                      |
|--------------------|----------------------|
| Species Reactivity | Human                |
| Class              | Primary              |
| Clonality          | Monoclonal           |
| Host               | Mouse                |
| Isotype            | Mouse igg1, κ        |
| Immunogen          | CD40 (BP50, TNFRSF5) |
| Clone              | G28.5                |
| Conjugate          | TRITC                |

**Biological Properties**

|             |                                                                                                      |
|-------------|------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                          |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | TRITC  |
| Excitation Wavelength | 544 nm |
| Emission Wavelength   | 570 nm |

**Applications**

G28.5 is an anti-human monoclonal antibody that targets the CD40 antigen. CD40 (also known as Bp50) is a 48 kD transmembrane protein that is expressed on the surface of cells like macrophages, dendritic cells and stem cells. CD40 is a component of key cellular pathways, in particular, the tumor necrosis factor-mediated signaling pathway, immune response-regulating cell surface receptor signaling pathway and CD40 signaling pathway. Also, in many organisms, it upregulates interleukin-12 production, is a positive regulator of isotype switching to IgG isotypes and is a promoter of protein phosphorylation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like CD40L, CD154 and TRAP. CD40 is a very popular antibody target, with over 30000 publications in the last decade.

CD40 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of neuroscience. This antibody was purified through affinity chromatography and conjugated to TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).