

TRITC Anti-human CD54 Antibody
R6-5-D6Catalog number: 105401I0, 105401I1
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 igg2a |
| Immunogen | CD54 (ICAM-1) |
| Clone | R6-5-D6 |
| 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

R6-5-D6 is an anti-human monoclonal antibody that targets the CD54 antigen. CD54 (also known as ICAM-1, ICAM1 or Ly-47) is a 85 - 110 kD member of the Ig superfamily that is expressed on the surface of cells such as B cells, endothelial cells and macrophages. CD54 is associated with a variety of biologically interesting macromolecules/ligands, for example, LFA-1, Mac-1 and Rhinovirus. CD54 is a fairly uncommon antibody target, with a little more than 4400 publications in the last decade. Even still, CD54 is essential for costimulatory molecules, neuroscience and immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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).