

iFluor™ 430 Anti-human CD54 Antibody
R6-5-D6Catalog number: 10540030, 10540031
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	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 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 iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSARIA™ III).