

**APC/iFluor™ 700 Anti-human/monkey  
CD154 Antibody \*5C8\***Catalog number: 115401F0, 115401F1, 115401F2  
Unit size: 25 tests, 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, monkey                    |
| Class              | Primary                          |
| Clonality          | Monoclonal                       |
| Host               | Mouse                            |
| Isotype            | Mouse igg2a                      |
| Immunogen          | CD154 (CD40L, gp39, TRAP, T-BAM) |
| Clone              | 5C8                              |
| Conjugate          | APC/iFluor™ 700                  |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                          |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | APC/iFluor™ 700 |
| Excitation Wavelength | 685 nm          |
| Emission Wavelength   | 710 nm          |

**Applications**

5C8 is an anti-human/monkey monoclonal antibody that forms an immune complex with the CD154 antigen. CD154 (also known as gp39) is a 32 - 39 kD member of the TNF superfamily that is expressed on the surface of cells such as T cells. CD154 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD40. CD154 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD154 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and

conjugated to APC/iFluor™ 700 (ex/em = 685/710 nm).