

**iFluor™ 670 Anti-human/monkey CD154
Antibody *5C8***Catalog number: 115400H0, 115400H1
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, monkey
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse igg2a
Immunogen	CD154 (CD40L, gp39, TRAP, T-BAM)
Clone	5C8
Conjugate	iFluor™ 670

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 670 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 670
Excitation Wavelength	671 nm
Emission Wavelength	682 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 iFluor™ 670 (ex/em = 671/682 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).