

**mFluor™ Red 700 Anti-mouse CD86  
Antibody \*GL-1\***Catalog number: 108600V0, 108600V1  
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 | Mouse            |
| Class              | Primary          |
| Clonality          | Monoclonal       |
| Host               | Rat              |
| Isotype            | Rat IgG2a kappa  |
| Immunogen          | CD86 (B7-2, B70) |
| Clone              | GL-1             |
| Conjugate          | mFluor™ Red 700  |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 700 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                    |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | mFluor™ Red 700 |
| Excitation Wavelength | 680 nm          |
| Emission Wavelength   | 695 nm          |

**Applications**

GL-1 is an anti-mouse monoclonal antibody that recognizes the CD86 antigen. CD86 (sometimes called B7-2 or B70) is a 80 kD glycoprotein that is expressed on the surface of cells such as T cells. CD86 is associated with a variety of biologically interesting macromolecules/ligands, namely, CD152 (CTLA-4). CD86 is a very popular antibody target, with over 21000 publications in the last decade. CD86 has been widely used in cell

biology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).