

**FITC Anti-mouse CD86 Antibody \*GL-1\***Catalog number: 108601I0, 108601I1  
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	FITC

**Biological Properties**

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

**Spectral Properties**

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 533/30 nm bandpass filter (for example, as in the BD FACSCanto™).