

**XFD647 Anti-mouse CD86 Antibody \*GL-1\***

Catalog Number: 10860180, 10860181

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	Lot specific (please consult certificate of analysis for given lot)
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	AF647

**Biological Properties**

---

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

**Spectral Properties**

---

Conjugate AF647

Excitation Wavelength 650 nm

Emission Wavelength 671 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 XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).