

XFD647 Anti-human CD85j Antibody *GHI/75*

Catalog Number: 10851180, 10851181

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2b kappa
Immunogen	CD85j (LILRB1, ILT2, LIR-1)
Clone	GHI/75
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

The GHI/75 monoclonal antibody recognizes human CD85j, a 110 kD member of the ILT/LIR family commonly found on the surface of natural killer cells and T cells. CD85j plays a role in important cellular pathways, in particular, the Fc receptor mediated inhibitory signaling pathway and immune response-inhibiting cell surface receptor signaling pathway. Moreover, it has been closely linked to critical biological processes like response to virus, especially defense response to virus. In some organisms, CD85j is a negative regulator of transforming growth factor-beta secretion, represses natural killer cell mediated cytotoxicity and inhibits dendritic cell apoptotic process. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands. CD85j is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD85j is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. 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).