

XFD647 Anti-human CD85 Antibody *17G10.2*

Catalog Number: 10850180, 10850181

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 IgG1 kappa
Immunogen	CD85g (LILRA4, ILT7)
Clone	17G10.2
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 17G10.2 monoclonal antibody binds to human CD85g, a single-pass type I membrane protein typically located on the surface of dendritic cells, neutrophils and eosinophils. CD85 plays a role in essential cellular pathways, in particular, the negative regulation of toll-like receptor 7 signaling pathway, negative regulation of toll-like receptor 9 signaling pathway and Fc-epsilon receptor signaling pathway. Moreover, in some organisms, it is a repressor of tumor necrosis factor production, is a negative regulator of toll-like receptor 9 signaling pathway and is involved in the negative regulation of interferon-alpha production. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands. CD85 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD85g is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of innate immunity and 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 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).