

PE/XFD700 Anti-human CD85 Antibody *17G10.2*

Catalog Number: 108501Q0,
108501Q1, 108501Q2
Unit Size: 25 tests, 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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
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	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).