

iFluor™ 430 Anti-human CD85 Antibody
17G10.2Catalog number: 10850030, 10850031
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD85g (LILRA4, ILT7)
Clone	17G10.2
Conjugate	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 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 iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).