

iFluor™ 594 Anti-human CD85j Antibody
GHI/75Catalog number: 108510C0, 108510C1
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 IgG2b kappa
Immunogen	CD85j (LILRB1, ILT2, LIR-1)
Clone	GHI/75
Conjugate	iFluor™ 594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 594
Excitation Wavelength	588 nm
Emission Wavelength	604 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 iFluor™ 594 (ex/em = 588/604 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).