

iFluor™ 488 Anti-human CD84 Antibody
CD84.1.21Catalog number: 10840050, 10840051
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 IgG2a kappa
Immunogen	CD84 (SLAMF5, Ly9b, GR6)
Clone	CD84.1.21
Conjugate	iFluor™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

CD84.1.21 is an anti-human monoclonal antibody that forms an immune complex with the CD84 antigen. CD84 (also known as Ly9b) is a 64 - 82 kD transmembrane glycoprotein that is found on the surface of cells like platelets, B cells, T cells, NK cells and dendritic cells. CD84 is associated with a variety of biologically interesting macromolecules/ligands, in particular, CD84. CD84 is a relatively rare antibody target, with fewer than

600 publications in the last decade. Even still, CD84 is typically 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™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 533/30 nm bandpass filter (for example, as in the BD FACSCanto™).