

Purified Anti-human CD84 Antibody
CD84.1.21

Catalog number: 10840000

Unit size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a kappa
Immunogen	CD84 (SLAMF5, Ly9b, GR6)
Clone	CD84.1.21

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.