

**mFluor™ Green 620 Anti-human CD148
Antibody *MEM-CD148/05***Catalog number: 114800U0, 114800U1
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
Immunogen	CD148 (DEP-1, p260)
Clone	MEM-CD148/05
Conjugate	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 nm

Applications

MEM-CD148/05 is an anti-human monoclonal antibody that forms an immune complex with the CD148 antigen. CD148 (sometimes called p260 or DEP-1) is a glycoprotein that is found on the surface of cells such as dendritic cells, platelets, B cells and NK cells. CD148 acts in key cellular pathways, for instance, the T cell receptor signaling pathway, negative regulation of epidermal growth factor receptor signaling pathway and

positive regulation of Fc-gamma receptor signaling pathway involved in phagocytosis. In addition, in many organisms, it upregulates MAPK cascade, is an inhibitor of cell growth and represses cell migration. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands. CD148 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD148 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 609/30 nm bandpass filter (for example, as in the Luminex Guava easyCyte).