

**PerCP Anti-rabbit/ cat/ human CD81
Antibody *M38***Catalog number: 108101V0, 108101V1, 108101V2
Unit size: 25 tests, 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	Rabbit, cat, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD81 (TAPA1, S5.7)
Clone	M38
Conjugate	PerCP

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 nm

Applications

The M38 monoclonal antibody binds with rabbit/ cat/ human CD81, a 26 kD member of the TM4SF tetraspanin family commonly found on the surface of nks, epithelial cells, T cells and natural killer cells. CD81 is associated with a variety of biologically interesting macromolecules/ligands, for example, CD21, leu-13 and CD19. CD81 is a fairly uncommon antibody target, with a little more than 6200 publications in the last decade. Even still, CD81 is vital to neuroscience and transcription factors research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 675/25 nm bandpass filter (for example, as in the BD Accuri™ C6 Plus).