

**iFluor™ 710 Anti-rabbit/ cat/ human CD81
Antibody *M38***Catalog number: 108100K0, 108100K1
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	Rabbit, cat, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD81 (TAPA1, S5.7)
Clone	M38
Conjugate	iFluor™ 710

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 710
Excitation Wavelength	717 nm
Emission Wavelength	739 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 iFluor™ 710 (ex/em = 717/739 nm).