

PerCP Anti-mouse CD105 Antibody
MJ7/18Catalog number: 110501S0, 110501S1, 110501S2
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD105 (Endoglin)
Clone	MJ7/18
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

MJ7/18 is an anti-mouse monoclonal antibody that recognizes the CD105 antigen. CD105 (sometimes called Endoglin) is a 90 kD glycoprotein that is expressed on the surface of cells like endothelial cells, stem cells and macrophages. CD105 is associated with a variety of biologically interesting macromolecules/ligands, in particular, TGF- β 3 and TGF- β 1. CD105 is a moderately popular antibody target, with over 11000 publications in the last decade. CD105 is essential for stem cells, immunology and angiogenesis research, frequently 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 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™

X-20).