

iFluor™ 860 Anti-mouse CD105 Antibody
MJ7/18Catalog number: 110500R0, 110500R1
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD105 (Endoglin)
Clone	MJ7/18
Conjugate	iFluor™ 860

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 860 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 860
Excitation Wavelength	853 nm
Emission Wavelength	878 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-β3 and TGF-β1. 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 iFluor™ 860 (ex/em = 853/878 nm).