

**iFluor™ 660 Anti-mouse CD105 Antibody**  
**\*MJ7/18\***Catalog number: 110500G0, 110500G1  
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™ 660

**Biological Properties**

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 660
Excitation Wavelength	663 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-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™ 660 (ex/em = 663/678 nm). It is compatible with the 640 nm laser and 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).