

XFD750 Anti-mouse CD105 Antibody *MJ7/18*

Catalog Number: 110501A0, 110501A1

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 | Lot specific (please consult certificate of analysis for given lot) |
| 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 | AF750 |

Biological Properties

| | |
|-----------------------|---|
| Appearance | Dark blue liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i> |

Spectral Properties

| | |
|-----------|-------|
| Conjugate | AF750 |
|-----------|-------|

Excitation Wavelength 752 nm

Emission Wavelength 776 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 XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).