

iFluor™ 700 Anti-human CD151 Antibody
50-6Catalog number: 115100J0, 115100J1
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD151 (PETA-3, SFA-1, GP27)
Clone	50-6
Conjugate	iFluor™ 700

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 700
Excitation Wavelength	690 nm
Emission Wavelength	713 nm

Applications

50-6 is an anti-human monoclonal antibody that is specific for the CD151 antigen. CD151 (alternatively called PETA-3, SFA-1 or GP27) is a 29 kD member of the tetraspanin family that is located on the surface of cells such as stem cells, epithelial cells, endothelial cells and platelets. CD151 is associated with a variety of biologically interesting macromolecules/ligands, namely, $\alpha 6$ and integrins $\alpha 3$. CD151 is a fairly uncommon

antibody target, with a little more than 1600 publications in the last decade. Even still, CD151 is vital to stem cells 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™ 700 (ex/em = 690/713 nm).