

## XFD700 Anti-human CD167a Antibody \*51D6\*

Catalog Number: 11670190, 11670191

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgM
Immunogen	CD167a (DDR1, trkE, CAK)
Clone	51D6
Conjugate	AF700

### Biological Properties

---

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 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	AF700
-----------	-------

Excitation Wavelength 696 nm

Emission Wavelength 719 nm

## Applications

---

51D6 is an anti-human monoclonal antibody that forms an immune complex with the CD167a antigen. CD167a (sometimes called DDR1) is a 5 kD glycoprotein that is found on the surface of cells such as dendritic cells and B cells. CD167a is associated with a variety of biologically interesting macromolecules/ligands, for example, FRS2, ShcA and collagens. CD167a is a relatively rare antibody target, with fewer than 10 publications in the last decade. Even still, CD167a is vital to immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD700 (ex/em = 696/719 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher).