

**iFluor™ 560 Anti-human CD167a Antibody**  
**\*51D6\***Catalog number: 116700A0, 116700A1  
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 IgM
Immunogen	CD167a (DDR1, trkE, CAK)
Clone	51D6
Conjugate	iFluor™ 560

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 560 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 560
Excitation Wavelength	560 nm
Emission Wavelength	571 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 iFluor™ 560 (ex/em =

560/571 nm). It is compatible with the 561 nm laser and 572/28 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).