

**mFluor™ Violet 510 Anti-human CD167a
Antibody *51D6***Catalog number: 11670110, 11670111
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	mFluor™ Violet 510

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 510 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 510
Excitation Wavelength	412 nm
Emission Wavelength	505 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 mFluor™ Violet 510 (ex/em = 412/505 nm). It is compatible with the 405 nm laser and 537/65 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).