

**mFluor™ Violet 450 Anti-human CD13
Antibody *WM15***Catalog number: 101300Z0, 101300Z1
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
Immunogen	CD13 (Aminopeptidase N, APN, gp150, ANPEP, PEPN)
Clone	WM15
Conjugate	mFluor™ Violet 450

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 nm

Applications

WM15 is an anti-human monoclonal antibody that targets the CD13 antigen. CD13 (sometimes referred to as gp150 or APN) is a 150 - 170 kD single-pass type II membrane protein that is found on the surface of cells such as epithelial cells, granulocytes, T cells, endothelial cells and macrophages. CD13 is associated with a variety of biologically interesting macromolecules/ligands, in particular, MEP1B, HNF1A, NGR and

Corona virus Receptor. CD13 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD13 has been widely used in stem cells and 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 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSCanto™ II).