

**mFluor™ Violet 450 Anti-human CD1  
Antibody \*SN13\***Catalog number: 100120Z0, 100120Z1  
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	CD1b (R1)
Clone	SN13
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**

The SN13 monoclonal antibody binds to human CD1b, a transmembrane protein commonly found on the surface of langerhans cells, dendritic cells and Tregs. In some organisms, CD1 is a positive regulator of T cell mediated cytotoxicity, and is associated with a variety of biologically interesting macromolecules/ligands, in particular,  $\beta$ -2-microglobulin. CD1 is a moderately popular antibody target, with over 15000 publications

in the last decade. CD1b is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. 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 457/45 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).