

**mFluor™ UV375 Anti-human CD138
Antibody *MI15***Catalog number: 113800X0, 113800X1
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	CD138 (Syndecan-1)
Clone	MI15
Conjugate	mFluor™ UV375

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ UV375 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ UV375
Excitation Wavelength	351 nm
Emission Wavelength	387 nm

Applications

The MI15 monoclonal antibody reacts with human CD138, a 100 - 200 kD transmembrane protein typically expressed on the surface of b cells, plasma cells, epithelial cells and endothelial cells. CD138 is involved with important cellular pathways, in particular, the cytokine-mediated signaling pathway and canonical Wnt signaling pathway. Also, in certain organisms, it enhances extracellular exosome assembly and is a

promoter of exosomal secretion. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands such as Fibronectin. CD138 is a fairly uncommon antibody target, with a little more than 7000 publications in the last decade. Even still, CD138 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of cell motility/cytoskeleton/structure, cell biology and synaptic biology. This antibody was purified through affinity chromatography and conjugated to mFluor™ UV375 (ex/em = 351/387 nm). It is compatible with the 355 nm laser and 379/28 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).