

iFluor™ 633 Anti-human CD138 Antibody
MI15Catalog number: 113800E0, 113800E1
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	iFluor™ 633

Biological Properties

Appearance	Black liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 633 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 633
Excitation Wavelength	640 nm
Emission Wavelength	654 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 iFluor™ 633 (ex/em = 640/654 nm). It is compatible with the 640 nm laser and 675/25 nm bandpass filter (for example, as in the BD Accuri™ C6 Plus).