

PE/XFD700 Anti-human CD138 Antibody *MI15*

Catalog Number: 113801P0,
113801P1, 113801P2
Unit Size: 25 tests, 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	Lot specific (please consult certificate of analysis for given lot)
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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).