

**mFluor™ Violet 510 Anti-human CD267
Antibody *1A1***Catalog number: 12670110, 12670111
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	N/a
Isotype	N/A
Immunogen	CD267 (TACI, TNFRSF13B)
Clone	1A1
Conjugate	mFluor™ Violet 510

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 510 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 510
Excitation Wavelength	412 nm
Emission Wavelength	505 nm

Applications

The 1A1 monoclonal antibody binds with human CD267, a 32 kD single-pass type iii membrane protein typically found on the surface of myeloma cells and B cells. CD267 is a component of important cellular pathways, namely, the tumor necrosis factor-mediated signaling pathway and cell surface receptor signaling pathway. Also, in some organisms, it is a suppressor of B cell proliferation. From a research standpoint, it is of

biological interest due to its association with important macromolecules/ligands like BAFF, BLYS and TALL1. CD267 is a relatively rare antibody target, with fewer than 50 publications in the last decade. Even still, CD267 has been widely used in costimulatory molecules and immunology research, frequently 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 510 (ex/em = 412/505 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant Analyzer 10).