

XFD488 Anti-human CD267 Antibody *1A1*

Catalog Number: 12670150, 12670151

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	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	N/a
Isotype	N/A
Immunogen	CD267 (TACI, TNFRSF13B)
Clone	1A1
Conjugate	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
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	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 527/32 nm bandpass filter (for example, as in the BD FACSMelody™).