

**iFluor™ 670 Anti-human CD267 Antibody**  
**\*1A1\***Catalog number: 126700H0, 126700H1  
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          | iFluor™ 670             |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 670 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 670 |
| Excitation Wavelength | 671 nm      |
| Emission Wavelength   | 682 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 iFluor™ 670 (ex/em = 671/682 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).