

**iFluor™ 405 Anti-human CD268 Antibody**  
**\*11C1\***Catalog number: 12680020, 12680021  
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 kappa         |
| Immunogen          | CD268 (BAFFR, TNFRSF13C) |
| Clone              | 11C1                     |
| Conjugate          | iFluor™ 405              |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Light yellow liquid                                                                                        |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 405 |
| Excitation Wavelength | 403 nm      |
| Emission Wavelength   | 427 nm      |

**Applications**

11C1 is an anti-human monoclonal antibody that forms an immune complex with the CD268 antigen. CD268 (alternatively called TNFRSF13C) is a 19 kD glycoprotein that is found on the surface of cells such as macrophages and granulocytes. CD268 is associated with a variety of biologically interesting macromolecules/ligands, for example, BAFF. CD268 is a relatively rare antibody target, with fewer than publications in

the last decade. Even still, CD268 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).