

**XFD594 Anti-human CD72 Antibody \*3F3,  
XFD594 Same Structure to Alexa Fluor™  
594\***Catalog number: 10720170, 10720171  
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 IgG2b                     |
| Immunogen          | CD72 ( Ly-19.2, Ly-32.2, Lyb-2) |
| Clone              | 3F3                             |
| Conjugate          | AF594                           |

**Biological Properties**

|             |  |
|-------------|--|
| Appearance  | Purple liquid  |
| Preparation | Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | AF594  |
| Excitation Wavelength | 590 nm |
| Emission Wavelength   | 618 nm |

**Applications**

The 3F3 monoclonal antibody binds with human CD72, a 39 - 43 kD transmembrane glycoprotein often located on the surface of B cells, dendritic cells, liver kupffer cells and macrophages. CD72 is associated with a variety of biologically interesting macromolecules/ligands, for

example, CD5. CD72 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD72 is commonly 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 594 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).