

**XFD680 Anti-human CD13 Antibody**  
**\*WM15, XFD680 Same Structure to Alexa**  
**Fluor™ 680\***

Catalog number: 10130190, 10130191  
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                                       |
| Immunogen          | CD13 (Aminopeptidase N, APN, gp150, ANPEP, PEPN) |
| Clone              | WM15   |
| Conjugate          | AF680  |

### Biological Properties

---

|             |  |
|-------------|--|
| Appearance  | liquid   |
| Preparation | Antibody purified by affinity chromatography and then conjugated with AF680 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

### Spectral Properties

---

|                       |        |
|-----------------------|--------|
| Conjugate             | AF680  |
| Excitation Wavelength | 681 nm |
| Emission Wavelength   | 704 nm |

### Applications

---

WM15 is an anti-human monoclonal antibody that targets the CD13 antigen. CD13 (sometimes referred to as gp150 or APN) is a 150 - 170 kD single-pass type II membrane protein that is found on the surface of cells such as epithelial cells, granulocytes, T cells, endothelial cells and

macrophages. CD13 is associated with a variety of biologically interesting macromolecules/ligands, in particular, MEP1B, HNF1A, NGR and Corona virus Receptor. CD13 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD13 has been widely used in stem cells and immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 680 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).