

## XFD488 Anti-human CD253 Antibody \*2E5\*

Catalog Number: 12530150, 12530151

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	Mouse
Isotype	Mouse IgG1
Immunogen	CD253 (TRAIL, Apo-2L, TNFSF10)
Clone	2E5
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

---

2E5 is an anti-human monoclonal antibody that recognizes the CD253 antigen. CD253 (sometimes called Apo-2L, TRAIL or TNFSF10) is a 30 kD member of the TNF ligand superfamily that is expressed on the surface of cells such as B cells and macrophages. CD253 is associated with a variety of biologically interesting macromolecules/ligands, for example, apo2. CD253 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD253 is essential for 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 525/50 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).