

iFluor™ 350 Anti-human CD253 Antibody
2E5Catalog number: 12530010, 12530011
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	CD253 (TRAIL, Apo-2L, TNFSF10)
Clone	2E5
Conjugate	iFluor™ 350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 350
Excitation Wavelength	345 nm
Emission Wavelength	450 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 iFluor™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).