

PerCP Anti-human CD253 Antibody *2E5*Catalog number: 125301V0, 125301V1, 125301V2
Unit size: 25 tests, 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	PerCP

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACSAria™ II).