

**PacBlue Anti-human CD340 Antibody**  
**\*24D2\***Catalog number: 134001K0, 134001K1  
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	CD340 (ERBB2, HER-2)
Clone	24D2
Conjugate	PacBlue

**Biological Properties**

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

**Spectral Properties**

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 nm

**Applications**

24D2 is an anti-human monoclonal antibody that targets the CD340 antigen. CD340 (alternatively called ERBB2 or HER-2) is a 185 kD transmembrane glycoprotein that is found on the surface of cells such as epithelial cells. CD340 has been closely linked to key biological processes like signal transduction, especially intracellular signal transduction. Furthermore, in certain organisms, it is an enhancer of cell growth,

is an enhancer of GTPase activity and plays a role in the upregulation of translation. CD340 is a member of important cellular pathways, for example, the cell surface receptor signaling pathway, transmembrane receptor protein tyrosine kinase signaling pathway and negative regulation of ERBB signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as EGFR. CD340 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD340 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of synaptic biology and immunology. This antibody was purified through affinity chromatography and conjugated to PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 445/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).