

PE/iFluor™ 750 Anti-human CD326 Antibody
323/A3Catalog number: 132601T0, 132601T1, 132601T2
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	CD326 (Ep-CAM)
Clone	323/A3
Conjugate	PE/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

The 323/A3 monoclonal antibody reacts with human CD326, a 40 kD glycoprotein commonly located on the surface of lungs and colons. In certain organisms, CD326 upregulates cell population proliferation, represses cell-cell adhesion mediated by cadherin and is an enhancer of transcription by RNA polymerase II. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands. CD326 is a fairly uncommon antibody target, with a little more than 1500 publications in the last decade. Even still, CD326 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell

biology and immunology. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 750 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 773/56 nm bandpass filter (for example, as in the Luminex Amnis CellStream).