

**iFluor™ 430 Anti-human CD324 Antibody
*67A4***Catalog number: 13240030, 13240031
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	CD324 (E-Cadherin)
Clone	67A4
Conjugate	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 nm

Applications

67A4 is an anti-human monoclonal antibody that targets the CD324 antigen. CD324 (also known as E-Cadherin) is a 100 kD member of the cadherin superfamily that is found on the surface of cells like erythrocytes. In certain organisms, CD324 is involved in the positive regulation of transcription, DNA-templated, enhances protein import into nucleus and negatively regulates cell migration. Also, it has been thought to be

involved with key biological processes such as cell-cell adhesion, specifically cell-cell adhesion mediated by cadherin. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands. CD324 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD324 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).