

**mFluor™ Blue 570 Anti-human CD24
Antibody *HI45***Catalog number: 102400T0, 102400T1
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	CD24 (Nectadrin, HSA)
Clone	HI45
Conjugate	mFluor™ Blue 570

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Blue 570
Excitation Wavelength	553 nm
Emission Wavelength	565 nm

Applications

HI45 is an anti-human monoclonal antibody that is specific for the CD24 antigen. CD24 (also known as Nectadrin or Ly-52) is a 35 - 45 kD transmembrane protein that is expressed on the surface of cells such as granulocytes, B cells and epithelial cells. CD24 acts in important cellular pathways, for instance, the immune response-regulating cell surface receptor signaling pathway, Wnt signaling pathway and intrinsic apoptotic

signaling pathway. Also, in some organisms, it is involved in the negative regulation of transforming growth factor beta3 production, plays a role in the upregulation of activated T cell proliferation and is involved in the positive regulation of MAP kinase activity. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like CD62P (P-Selectin). CD24 is a moderately popular antibody target, with over 12000 publications in the last decade. CD24 has a variety of applications in cell biology and immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).