

iFluor™ 820 Anti-human CD70 Antibody
Ki-24Catalog number: 107000P0, 107000P1
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 IgG3 kappa
Immunogen	CD70 (CD27 ligand, CD27L, TNFSF7)
Clone	Ki-24
Conjugate	iFluor™ 820

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 820 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 nm

Applications

The Ki-24 monoclonal antibody binds to human CD70, a 50 kD member of the tumor necrosis factor superfamily frequently found on the surface of T cells, dendritic cells, B cells and natural killer cells. CD70 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD27 ligand. CD70 is a fairly uncommon antibody target, with a little more than 2900 publications in the last decade. Even still, CD70

has been widely used in immunology and costimulatory molecules research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 820 (ex/em = 822/850 nm).