

**mFluor™ Blue 570 Anti-human CD117
Antibody *104D2***Catalog number: 111700T0, 111700T1
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	CD117 (c-kit, SCFR)
Clone	104D2
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

The 104D2 monoclonal antibody reacts with human CD117, a 145 kD single-pass type i membrane protein typically found on the surface of leukemias, embryonic stem cells, mesenchymal stem cells and mast cells. CD117 has been closely linked to vital biological processes like germ cell migration, specifically ectopic germ cell programmed cell death. In addition, in some organisms, it acts to positively regulate pseudopodium

assembly, plays a role in the upregulation of long-term neuronal synaptic plasticity and is an inhibitor of programmed cell death. CD117 is a component of critical cellular pathways, for example, the transmembrane receptor protein tyrosine kinase signaling pathway, Kit signaling pathway and Fc receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as MGF, SCF and KL. CD117 is a moderately popular antibody target, with over 11000 publications in the last decade. CD117 has been widely used in stem cells and immunology research, frequently 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 582/15 nm bandpass filter (for example, as in the BD FACSaria™ III).