



# CD117-ECD

# PN B38307

## Antibody

Specificity: Human  
Clone: 104D2D1  
Isotype: IgG1 kappa mouse

## Reagent

Line: IOTest  
Format: Liquid  
Size: 0.5 mL  
Status: ASR

## Antibody characteristics

The monoclonal antibody 104D2D1 was assigned to CD117 during the 6th HLDA Workshop on Human Leucocyte Differentiation Antigens held in Kobe, Japan, in 1996 (WS Code: C-30. Section C).

## Antigen characteristics

Other name: c-kit, stem cell factor receptor (SCFR)      Gene Name: KIT  
Gene Location: 4q11-q12      Molecular Weight: 145 kDa

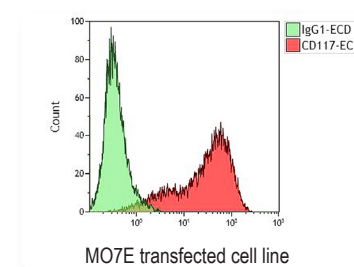
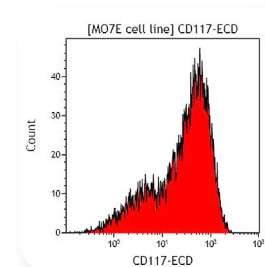
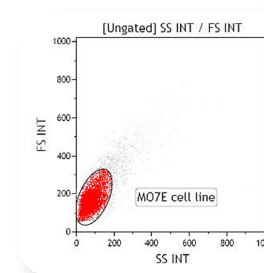
The CD117 antigen, also known as Stem Cell Factor Receptor (SCFR), mast-cell-Kit, and steel factor receptor, is a transmembrane glycoprotein encoded by the c-kit proto-oncogen. The CD117 molecule belongs to the class III Receptor Tyrosine Kinase (RTK) family. Within the hematopoietic compartment, the CD117 molecule is expressed on approximately 50% of CD34<sup>+</sup> progenitors engaged in erythrocytic, myelomonocytic and megakaryocytic differentiation. Although CD117 is primarily a marker for non-lymphoid progenitor, it has been reported to be detected on early lymphoid progenitor. CD117 expression is found on a small subset of resting NK cells (CD56 bright), and about 30% of immature CD3<sup>-</sup> CD4<sup>-</sup> CD8<sup>-</sup> thymocytes. CD117 is also expressed on mast cells and detected on non-hematopoietic cells such as reproductive system, melanocytes and embryonic brain.

## Formats available

	Volume	Line	Status	Part #		Volume	Line	Status	Part #
PE	2 mL	IOTest	ASR	IM2732U	PC7	1 mL	IOTest	ASR	IM3698
ECD	0.5 mL	IOTest	ASR	B38307	APC	1 mL	IOTest	ASR	IM3638
PC5	1 mL	IOTest	ASR	IM2733U	APC-A750 <sup>(1)</sup>	0.5 mL	IOTest	ASR	A86051
PC5.5	0.5 mL	IOTest	ASR	A66333					

(1) APC-Alexa Fluor\* 750

## Example of results



MO7E transfected cell line

## References

- ◆ Uoshima, N., Ozawa, M., Kimura, S., Tanaka, K., Wada, K., Kobayashi, Y., Kondo, M., «Changes in c-Kit expression and effects of SCF during differentiation of human erythroid progenitor cells», 1995, Br. J. Haematol., 91, 30-36.
- ◆ Ashman, L.K., Cambareli, A.C., Nguyen, L., Bühring, H.-J., «CD117 Workshop panel report», 1997, Leucocyte Typing VI, White Cell Differentiation Antigens. Kishimoto, T., et al, Eds., Garland Publishing, Inc., 816-818.
- ◆ Kikutani, H., Kishimoto, T., «The cytokine receptors: Section report», 1995, Leucocyte Typing V, White Cell Differentiation Antigens. Schlossman, S.F., et al., Eds., Oxford University Press, 1855-1864.

ASR: Analyte Specific Reagent. Analytical and performance characteristics are not established.

\*Alexa Fluor is registered trademark of Molecular Probes, Inc.  
Beckman Coulter, the stylized logo and IOTest are trademarks of Beckman Coulter, Inc. and are registered in the USPTO.