

**ANTI-HUMAN CD117 PE**

**PE-Conjugated Mouse Monoclonal Antibody Against Human CD117  
Clone 104D2**

**Catalog # 10531**

**50 tests**

**SPECIFICITY:**

CD117 is the receptor for stem cell factor (SCF), c-Kit. CD117 is expressed on ~1-4% of bone marrow cells, including hematopoietic stem cells, pluripotent, myelomonocytic and erythroid progenitor cells and erythroblasts, CD117 is also expressed on tissue mast cells, melanocytes, primordial germ cells and acute myeloid leukemia cells. Anti-CD117 antibody clone 104D2 does not block the epitope that binds SCF.

**CLONE:** 104D2

**ISOTYPE:** IgG<sub>1</sub>κ (mouse)

**FORMAT:**

PE-conjugated antibody in 1 mL Phosphate Buffered Saline (PBS) containing gelatin and 0.1% (w/v) sodium azide.

**STABILITY AND STORAGE:**

Store at 4°C. Do not freeze. Product is stable for at least 6 months.

**APPLICATIONS AND DIRECTIONS FOR USE:**

**Flow cytometry:**

Recommended amount per 1x10<sup>6</sup> cells in a volume of 100 µL: 20 µL  
Appropriate conditions should be established for each application.

**Cell separation:**

Positive selection of CD117 (Kit)<sup>+</sup> cells with StemCell's reagents for immunomagnetic cell separation. Please contact us for more information.

**NOT FOR CLINICAL USE; INCLUDING IN VITRO DIAGNOSTIC USE, AND EX VIVO OR IN VIVO THERAPEUTIC USE IN CLINICAL TRIALS OR IN CLINICAL PRACTISE.  
THIS REAGENT IS FOR RESEARCH ONLY.**

***Hazardous Ingredient: Sodium Azide.*** Avoid exposure to skin and eyes, ingestion and contact with heat, acids and metals. Wash exposed skin with soap and water. Flush eyes with water. Dilute with running water before discharging into plumbing.

**REFERENCES:**

1. Barclay AN, Brown MH, Law SKA, McKnight AJ, Tomlinson MG, van der Merwe PA, eds. 1997. The Leukocyte Antigens Facts Book, 2<sup>nd</sup> Edition, CD117. Academic Press, NY, p 394-396.
2. Kishimoto T et al. Eds.1998. Leukocyte Typing VI. White Cell Differentiation Antigens. Garland Publishing Inc, New York NY, p1191-1192.
3. Olweus J, Terstappen LW, Thompson PA, Lund-Johansen F. Expression and function of receptors for stem cell factor and erythropoietin during lineage commitment of human hematopoietic progenitor cells. Blood 88: 1594-607, 1996.
4. Wognum AW, de Jong MO, Wagemaker G. Differential expression of receptors for hemopoietic growth factors on subsets of CD34+ hemopoietic cells. Leuk Lymphoma 24:11-25, 1996.