

**ANTI-HUMAN CD19**

**Conjugated Mouse Monoclonal Antibodies Against Human CD19  
Clone 4G7**

<b>FITC Conjugate:</b>	<b>Catalog # 10409</b>	<b>100 tests</b>
<b>PE Conjugate:</b>	<b>Catalog # 10509</b>	<b>50 tests</b>

**SPECIFICITY:**

CD19 is expressed on all stages of B-cell differentiation (from CD34<sup>+</sup> pre B-cells to mature B-cell), but is lost upon maturation to plasma cells. CD19 is also expressed on follicular dendritic cells.

**CLONE:** 4G7

**ISOTYPE:** IgG<sub>1</sub>κ (mouse)

**FORMAT:**

FITC-conjugate: 50 µg in 2 mL (25 µg/mL) in phosphate buffered saline (PBS) containing gelatin and 0.1% (w/v) sodium azide.

PE-conjugate : 25 µg in 1 mL (25 µg/mL) in 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:

- FITC-conjugate (Cat No. 10409): 20 µL
- PE-conjugate (Cat No. 10509): 20 µL

Appropriate conditions should be established for each application.

**Cell separation:**

Positive selection of CD19<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, CD19. Academic Press, NY, p 179-180.
2. Kishimoto T et al. Eds.1998. Leukocyte Typing VI. White Cell Differentiation Antigens. Garland Publishing Inc, New York NY, p1123-1124.