

ANTI-MOUSE CD4

**Conjugated Rat Monoclonal Antibodies Against Mouse CD4
Clone H129.19**

FITC-conjugate:	Catalog # 10701	100 mg
	Catalog # 10702	500 mg
PE-conjugate:	Catalog # 10801	100 mg
	Catalog # 10802	200 mg

SPECIFICITY:

CD4/L3T4 is a 52 kDa differentiation antigen expressed on most thymocytes and a subpopulation of mature T-lymphocytes, i.e., the helper/inducer subpopulation. In the mouse, CD4 is not expressed on monocytes and macrophages. The H129.19 antibody reacts with all mouse strains tested. The H129.19 monoclonal antibody blocks binding of the anti-mouse CD4 monoclonal antibody GK1.5 and inhibits T helper cell responses to antigenic or mitogenic stimuli.

CLONE: H129.19

ISOTYPE: IgG_{2a}K (rat)

FORMAT:

FITC-conjugate: Catalog # 10701: 100 µg in 0.2 mL (0.5 mg/mL) diluent
Catalog # 10702: 500 µg in 1.0 mL (0.5 mg/mL) diluent
PE-conjugate: Catalog # 10801: 100 µg in 0.5 mL (0.2 mg/mL) diluent
Catalog # 10802: 200 µg in 1.0 mL (0.2 mg/mL) diluent
Diluent: Aqueous buffered solution, containing 0.09% (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 1x10⁶ cells in a volume of 100 µL:

- FITC-conjugate (Cat No. 10701,10702): ≤ 1 µg (2 µL)
- PE-conjugate (Cat No. 10801, 10802): ≤ 1 µg (5 µL)

Appropriate conditions should be established for each application.

Cell separation:

Positive selection of CD4+ 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 ON**

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, 2nd Edition, CD4. Academic Press, NY, p 141-142.
2. Dialynas DP, Quan ZS, Wall KA, Pierres A, Quintans J, Loken MR, Pierres M, Fitch FW. Characterization of the murine T cell surface molecule, designated L3T4, identified by monoclonal antibody GK1.5: similarity of L3T4 to the human Leu-3/T4 molecule. 1983, J Immunol 13:2445-2451.
3. Saizawa K, Haque S, Jones B, Rojo J, Tite JP, Kaye J, Janeway CA Jr . The L3T4 molecule is part of the helper T-cell antigen/Ia recognition complex. 1987, Ann Inst Pasteur Immunol 138: 138-143