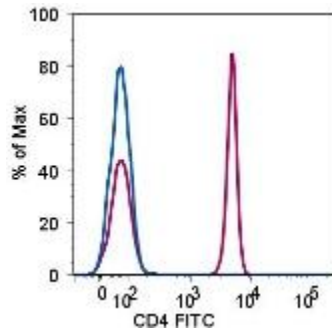


## CD4 FITC

**Catalog Number:** 8011-0047

Also known as: Leu-3a

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



Fluorescence profiles of normal human peripheral blood lymphocytes unstained (blue histogram) or stained with CD4 (Leu-3a) FITC (purple histogram).

### Product Information



**Contents:** CD4 FITC

**Catalog Number:** 8011-0047

**Clone:** SK3

**Concentration:** 5 uL (0.06 ug)/test (a test is defined as the amount that will stain 1 x 10<sup>6</sup> cells in 100 uL)

**Host/Isotype:** Mouse IgG1, kappa

**HLDA Workshop:** I

**Formulation:** aqueous buffer, 0.09% sodium azide, may contain carrier protein/stabilizer



**Temperature Limitation:** Store at 2-8°C. Do not freeze. Light-sensitive material.



**Batch Code:** Refer to vial



**Analyte Specific Reagent**



**Manufacturer:** eBioscience, Inc., 10255 Science Center Drive, San Diego, CA 92121, USA



**Use By:** Refer to vial

**Caution, contains Azide**

### Description

The SK3 monoclonal antibody reacts with human CD4, a 59-kDa cell surface receptor expressed by a majority of thymocytes, a subpopulation of mature T helper cells, and, in low levels, monocytes. CD4 is a receptor for the human immunodeficiency virus (HIV).

### References

Knapp W, Dorken B, et al, eds. Leucocyte Typing IV: White Cell Differentiation Antigens. Oxford University Press. New York. 1989.

Bernard A, Boumsell L, Hill C. Joint report of the first international workshop on human leucocyte differentiation antigens by the investigators of the participating laboratories. In: Bernard A, Boumsell L, Dausset J, Milstein C, Schlossman SF, eds. Leucocyte Typing. New York, NY: Springer-Verlag; 1984:9-108.