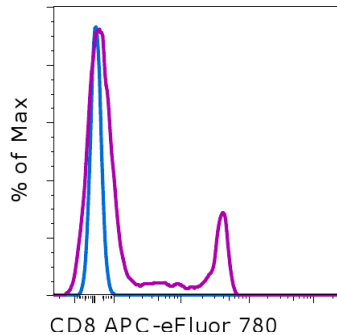


## CD8 APC-eFluor® 780

**Catalog Number:** 8047-0087

Also known as: Leu-2a

**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 CD8 APC-eFluor® 780 (purple histogram).

### Product Information



**Contents:** CD8 APC-eFluor® 780

**Catalog Number:** 8047-0087

**Clone:** SK1

**Concentration:** 5 uL (0.5 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 SK1 monoclonal antibody reacts with the human CD8 alpha molecule, an approximately 32-34 kDa cell surface receptor expressed either as a heterodimer with the CD8 beta chain (CD8 alpha/beta) or as a homodimer (CD8 alpha/alpha). A majority of thymocytes and a subpopulation of mature T cells and NK cells express CD8. CD8 binds to MHC class I and through its association with protein tyrosine kinase p56 lck plays a role in T-cell development and activation of mature T cells.

### References

Campanelli R, Palermo B, et al. Human CD8 co-receptor is strictly involved in MHC-peptide tetramer-TCR binding and T cell activation. *International Immunology*. 2002 May; 14(1):39-44.

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.