

Validation data for Anti-mCD8-mIgG2a

<https://www.invivogen.com/anti-mcd8-migg2a-invivoft>

For research use only

Version 20K10-NJ

Anti-mCD8-mIgG2a InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) designed for *in vivo* depletion in mice. This mAb features the variable region of the previously described anti-mCD8 (YTS 169.4 clone, rat IgG2b κ) mAb and a murine IgG2a constant region. The binding capacity of Anti-mCD8-mIgG2a InvivoFit™ to the murine CD8 (mCD8) antigen has been confirmed using cell surface staining with EL4 cells stably expressing mCD8 (EL4-mCD8) (Figure 1). The *in vivo* depletion of peripheral blood CD8⁺ T cells using Anti-mCD8-mIgG2a InvivoFit™ has been confirmed (Figure 2).

Validation of Anti-mCD8-mIgG2a binding by FACS

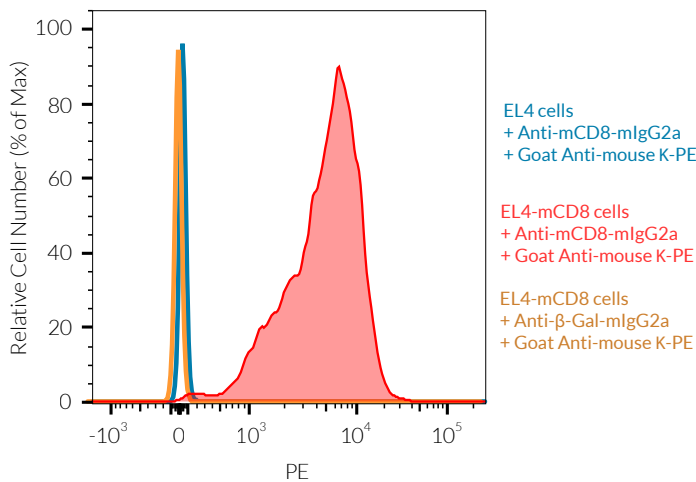


Figure 1: Cell surface staining of murine CD8 using Anti-mCD8-mIgG2a InvivoFit™ mAb.

~5x10⁵ EL4 (parental) or EL4-mCD8 cells were incubated with 2 μ g of Anti-mCD8-mIgG2a InvivoFit™ mAb or an isotype control for 1h at 4°C. Cells were then washed and incubated with 0.25 μ g of goat anti-mouse κ -light chain coupled to PE for 1h at 4°C. Cell surface staining was analyzed by flow cytometry.

Validation of *in vivo* CD8⁺ T cell depletion using Anti-mCD8-mIgG2a InvivoFit™

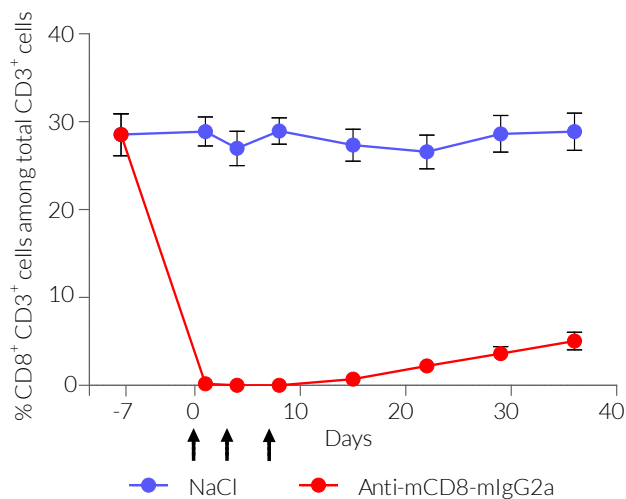


Figure 2: Analysis of CD8⁺ T cell population in mice peripheral blood upon injection of Anti-mCD8-mIgG2a InvivoFit™ mAb.

Blood samples (20 μ l) from SWR/J mice aged 10 weeks were harvested 1 week prior to a series of intraperitoneal (IP) injections with the depleting Anti-mCD8-mIgG2a InvivoFit™ mAb (200 μ g) or NaCl (control), and 24h after each IP injection. IP injections were performed at day 0, +3, and +7 (arrows). Thereafter, blood samples were harvested once a week. To assess the kinetics of the *in vivo* depletion of CD8⁺ T cells, peripheral blood was stained with the following conjugated antibodies: Anti-CD45-Pe-Cy7 (clone 30-F11), Anti-CD3e-PE (clone 145-2C11), Anti-CD4-FITC (clone RMA-4), and Anti-CD8-APC (clone H35-17.2) according to standard procedures. Cell surface staining was analyzed by flow cytometry. Data are represented as the frequency of CD8⁺ T cells among the total CD3⁺ T cell population.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3622-3480
E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com