

Validation data for Anti-mTIGIT-mIgG2a

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

For research use only

Version 21B01-NJ

Anti-mTIGIT-mIgG2a InvivoFit™ is an Anti-mTIGIT mAb featuring the variable region of the previously described 10A7 hamster IgG2a clone and a murinized IgG2a constant region. The cell surface staining capacity of Anti-mTIGIT-mIgG2a InvivoFit™ has been confirmed with EL4 cells stably expressing mTIGIT (EL4-mTIGIT) (Figure 1).

Validation of Anti-mTIGIT-mIgG2a InvivoFit™ by FACS

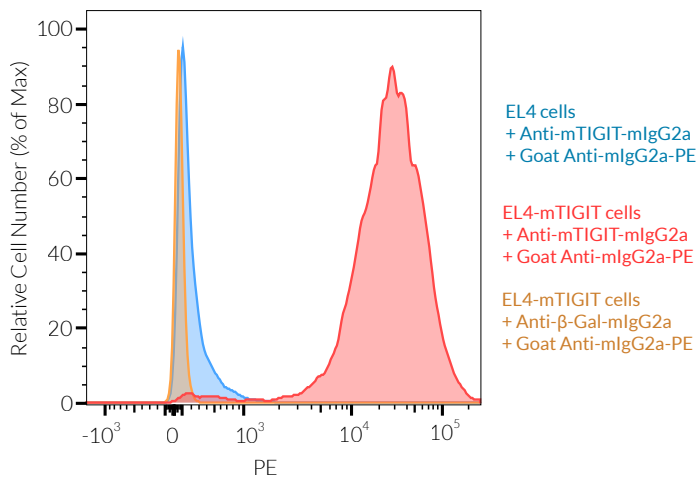


Figure 1: Cell surface staining of murine TIGIT using Anti-mTIGIT-mIgG2a InvivoFit™ mAb.

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

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