

# Validation data for Anti-mCTLA4-mIgG2a

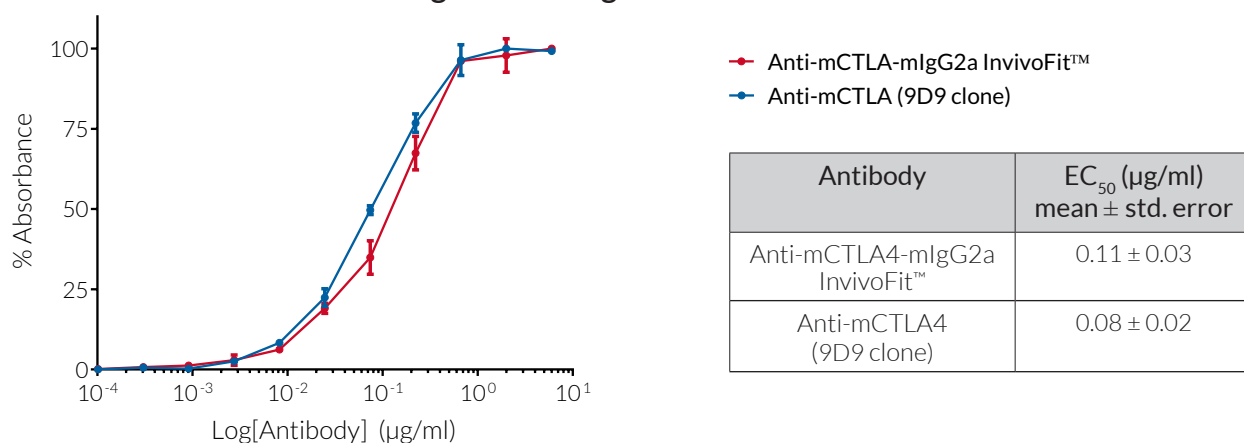
<https://www.invivogen.com/anti-mctla4-migg2a-invivofit>

For research use only

Version 19D23-MM

Anti-mCTLA4-mIgG2a InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) designed for *in vivo* studies in mice. This mAb features the variable region of the previously described anti-mCTLA (9D9 clone) mAb and a murine IgG2a constant region. The anti-mCTLA (9D9 clone) mAb recognizes murine CTLA-4 (mCTLA-4). It blocks CTLA-4-mediated negative signals that physiologically downregulate T cell activation, and thereby unleashes T cell activity. The binding capacity of Anti-mCTLA4-mIgG2a InvivoFit™ to the mCTLA-4 antigen has been confirmed using ELISA.

## Evaluation of Anti-mCTLA4-mIgG2a binding to mCTLA-4



ELISA binding of Anti-mCTLA4-mIgG2a InvivoFit™ to recombinant murine CTLA-4. A dilution series of Anti-mCTLA4-mIgG2a InvivoFit™ and anti-mCTLA4 (9D9 clone) was used for the detection of coated mCTLA-4 antigen at 5 µg/ml. Detection of bound antibody was performed using an HRP-conjugated anti-murine  $\kappa$ -light chain secondary antibody. HRP activity was assessed by a colorimetric assay using Tetramethylbenzidine (TMB) substrate. Optical density was measured at 490 nm. Data are shown as percentage absorbance.

### 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](mailto:info@invivogen.com)