

Validation data for Anti-mPD-1-mIgG1e3 InvivoFit™

<https://www.invivogen.com/anti-mpd1-migg1e3-invivofit>

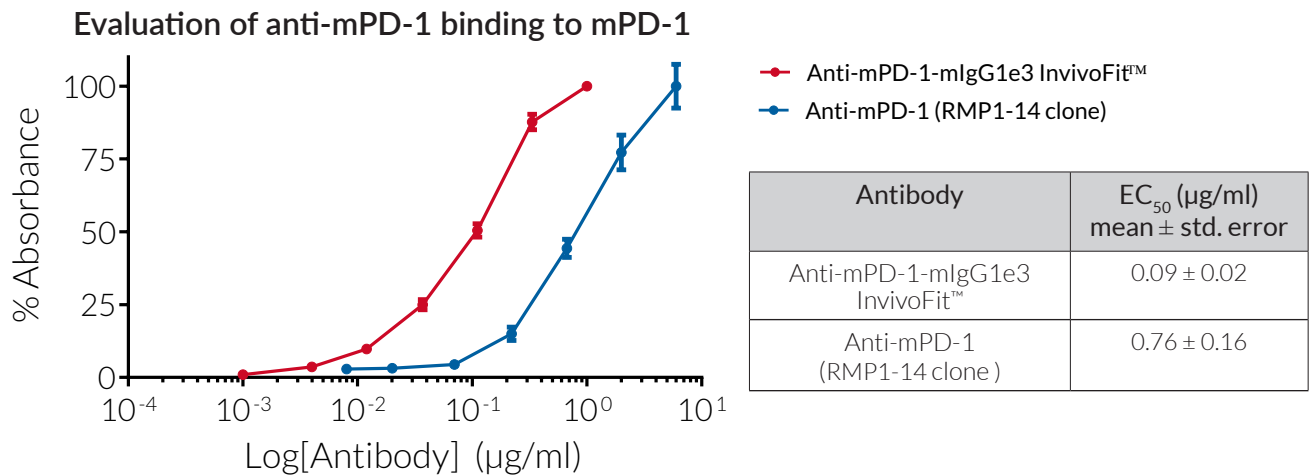
For research use only

Version 19D24-MM

Anti-mPD-1-mIgG1e3 InvivoFit™ is a recombinant monoclonal antibody (mAb) designed for *in vivo* studies in mice. This mAb features the variable region of the previously described anti-mPD-1 RMP1-14 mAb and the engineered murine IgG1e3 constant region. The RMP1-14 mAb targets and blocks murine programmed cell death 1 (mPD-1). However, the RMP1-14 mAb, along with other commercially available anti-mPD-1 mAbs, were generated in rats, and are therefore immunogenic in mice. Additionally, these mAbs feature the IgG2a isotype which promotes antibody-dependent cellular cytotoxicity.

Anti-mPD-1-mIgG1e3 InvivoFit™ was generated by recombinant DNA technology so that it is ~85% murine (constant region) and with a point mutation D265A (a replacement of aspartic acid by alanine at position 265), resulting in the complete loss of cytolytic effector function. Anti-mPD-1-mIgG1e3 InvivoFit™ is thus ideal for blocking the mPD-1 receptor without causing T cell depletion.

The binding capacity of Anti-mPD-1-mIgG1e3 InvivoFit™ to mPD-1 has been confirmed using ELISA.



ELISA binding of Anti-mPD-1-mIgG1e3 InvivoFit™ to recombinant murine PD-1. A dilution series of Anti-mPD-1-mIgG1e3 InvivoFit™ and anti-mPD-1 (RMP1-14 clone) was used for the detection of coated mPD-1 antigen at 2 µg/ml. Detection of bound antibody was performed using an HRP-conjugated anti-murine IgG secondary antibody for Anti-mPD-1-mIgG1e3 InvivoFit™ and an HRP-conjugated anti-rat IgG2a secondary antibody for anti-mPD-1 (RMP1-14 clone). 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

 **InvivoGen**
www.invivogen.com