

Validation data for Anti-mgp75-mlgG2a

<https://www.invivogen.com/anti-mgp75-migg2a-invivofit>

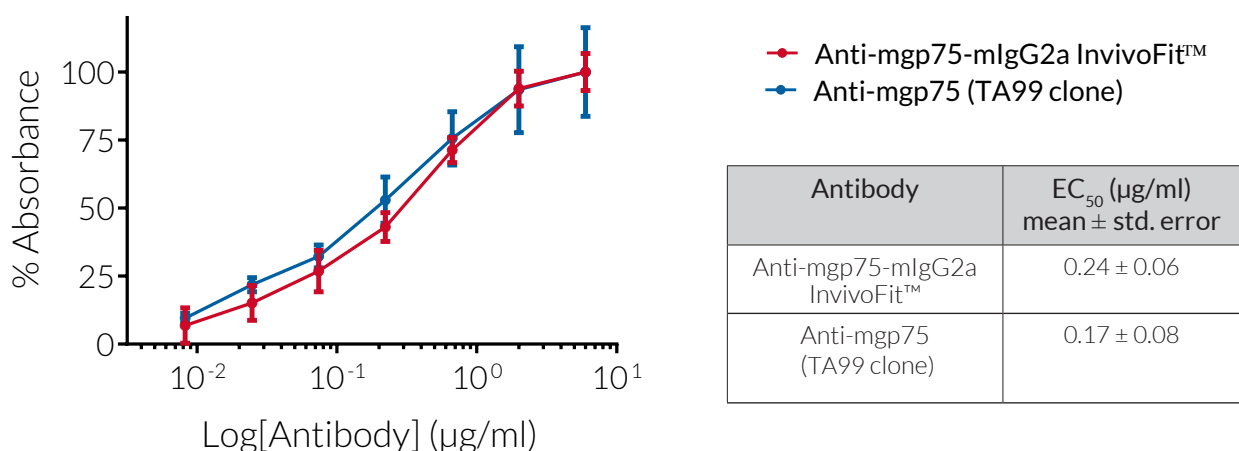
For research use only

Version 19C23-MM

Anti-mgp75-mlgG2a InvivoFit™ is a recombinant mouse anti-mouse monoclonal antibody (mAb) designed for *in vivo* studies in mice. This mAb features the variable region of the previously described anti-mgp75 hybridoma clone TA99 and the murine IgG2a constant region. The TA99 clone reacts with both human and mouse glycoprotein 75 (gp75), the mature form of tyrosinase-related protein 1 (TYRP1).

Anti-mgp75-mlgG2a InvivoFit™ has a sequence that is 100% murine (constant and variable regions), thus ensuring minimal immunogenicity upon repeated injections into mice. Notably, IgG2a is the isotype of choice for this antibody, as anti-mgp75 mAbs featuring the IgG2a constant regions display enhanced tumor clearance and antibody-dependent cellular cytotoxicity. The binding capacity of Anti-mgp75-mlgG2a InvivoFit™ to the murine gp75 (TYRP1) antigen has been confirmed using ELISA.

Evaluation of Anti-mgp75-mlgG2a binding to mgp75 (TYRP1)



ELISA binding of Anti-mgp75-mlgG2a InvivoFit™ to recombinant murine gp75 protein. A dilution series of Anti-mgp75-mlgG2a InvivoFit™ and anti-mgp75 (TA99 clone) was used for the detection of coated mgp75 (mTYRP1) antigen at 2 µg/ml. Detection of bound antibody was performed using an HRP-conjugated anti-murine IgG2a 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