

Validation data for Anti-mGITR-mIgG2a InvivoFit™

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

For research use only

Version 23C20-AK

Anti-mGITR-mIgG2a InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) featuring the variable region of the previously described anti-mGITR clone DTA-1. GITR (Glucocorticoid-Induced TNFR-Related protein) is a co-stimulatory immune checkpoint protein found on a variety of immune cells, and is highly involved in the pathogenesis of tumor, inflammation and autoimmune diseases. The ability of Anti-mGITR-mIgG2a InvivoFit™ to bind the murine GITR has been validated by ELISA, using coated mGITR and an anti-mIgG-HRP secondary antibody (Figure 1).

Validation of Anti-mGITR-mIgG2a InvivoFit™ by ELISA

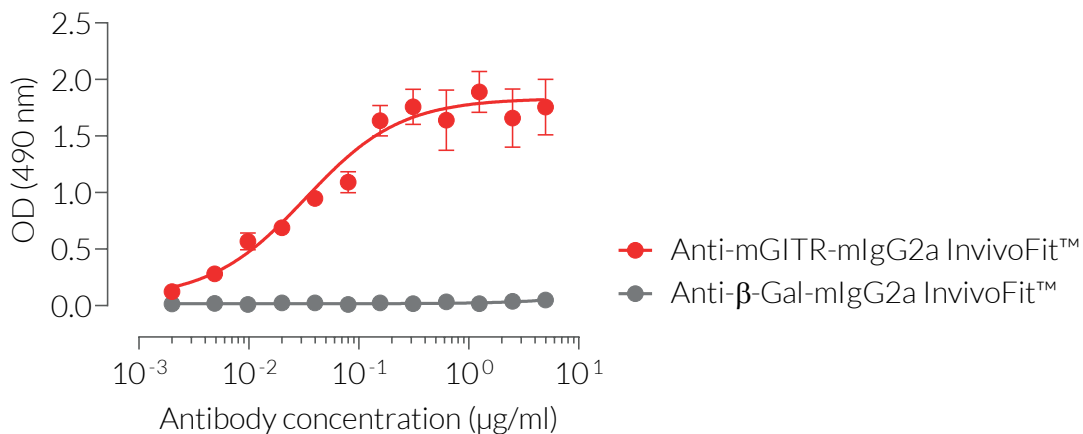


Figure 1. Binding of Anti-mGITR-mIgG2a InvivoFit™ mAb to coated mouse GITR protein.

Murine GITR (5 µg/ml) was coated on ELISA plates overnight. A 2-fold serial dilution of the Anti-mGITR-mIgG2a InvivoFit™ (red curve) or of the Anti-β-Gal-mIgG2a control antibody (grey curve) was performed for the capture step. An HRP-labelled anti-mIgG antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Data are shown as optical density (OD) at 490 nm (mean ± SEM).

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 Asia: +852 3622-3480
E-mail: info@invivogen.com