

# Validation data for Anti-mIL-16-mIgG1e3 InvivoFit™

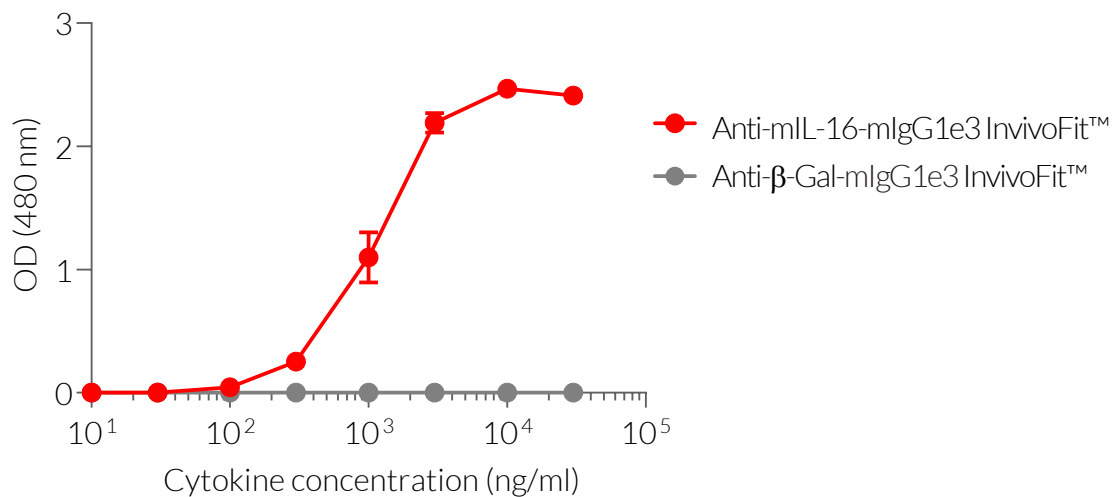
<https://www.invivogen.com/anti-mil16-igg1e3>

For research use only

Version 23E09-AK

Anti-mIL-16-mIgG1e3 InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) designed to detect murine interleukin 16 (mIL-16). This pleiotropic cytokine plays an important role in modulating T cell activation, viral defense, and inflammatory diseases. The ability of Anti-mIL-16-mIgG1e3 InvivoFit™ to bind the murine IL-16 has been validated by ELISA, using coated mIL-16 and an anti-mIgG-HRP secondary antibody (Figure 1).

## Validation of Anti-mIL-16-mIgG1e3 InvivoFit™ by ELISA



**Figure 1. Binding of Anti-mIL-16-mIgG1e3 InvivoFit™ mAb to coated mIL-16 protein.** Coated ELISA plates with increasing concentrations of mIL-16 were incubated with 2 µg/ml of Anti-mIL-16-mIgG1e3 InvivoFit™ mAb (red curve) or Anti-β-Gal-mIgG1e3 InvivoFit™ control antibody (grey curve) for 2h, followed by an HRP-conjugated anti-mouse secondary antibody (1/1000 dilution). Data are shown as optical density (OD) at 480 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](mailto:info@invivogen.com)