

Validation data for Anti-mIL-16-mIgG1e3

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

For research use only

Version 23E09-AK

Anti-mIL-16-mIgG1e3 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 to bind the mIL-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 by ELISA

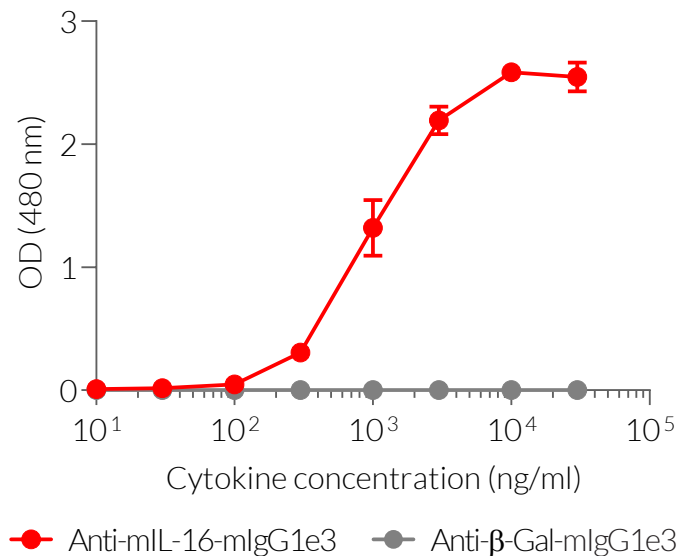


Figure 1. Binding of Anti-mIL-16-mIgG1e3 mAb to coated mIL-16 protein. Coated ELISA plates with increasing concentrations of mIL-16 were incubated with 2 µg/ml of Anti-mIL-16-mIgG1e3 mAb (red curve) or Anti-β-Gal-mIgG1e3 control antibody (grey curve) for 2h, followed by an HRP-conjugated anti-mouse secondary 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