

Validation data for Anti-hIL-16-hIgG1

<https://www.invivogen.com/anti-hIL16-igg>

For research use only

Version 23E09-AK

Anti-hIL-16-hIgG1 is a recombinant monoclonal antibody (mAb) designed to detect human interleukin 16 (hIL-16). This pleiotropic cytokine plays an important role in modulating T cell activation, viral defense, and inflammatory diseases. The ability of Anti-hIL-16-hIgG1 to bind hIL-16 has been validated by ELISA (Figure 1).

Validation of Anti-hIL-16-hIgG1 by ELISA

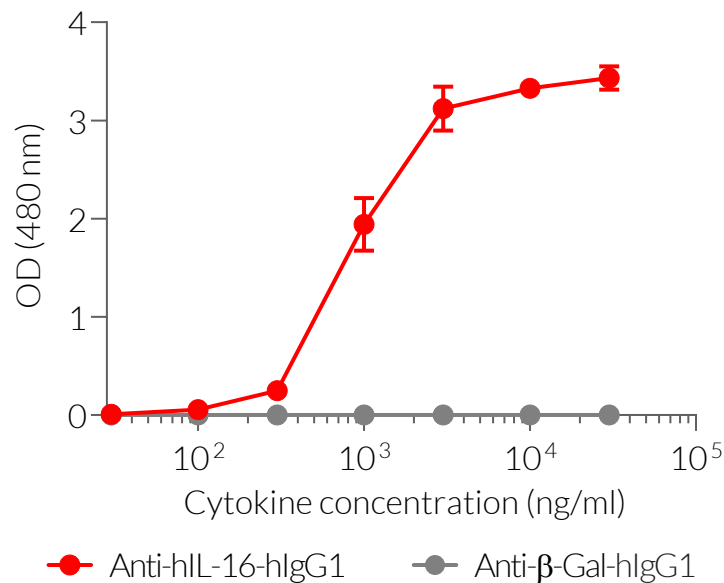


Figure 1. Binding of Anti-hIL-16-hIgG1 mAb to coated hIL-16 protein. Coated ELISA plates with increasing concentrations of hIL-16 were incubated with 2 µg/ml of Anti-hIL-16-hIgG1 mAb (red curve) or Anti-β-Gal-hIgG1 control antibody (grey curve) for 2h, followed by an HRP-conjugated anti-human secondary 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