

# Validation data for recombinant human IL-16

<https://www.invivogen.com/human-mouse-il16>

For research use only

Version 23E16-NJ

Interleukin 16 (IL-16) is a pro-inflammatory cytokine and alarmin. Its activity is mediated by the release of its C-terminal portion, after cleavage by caspase-3. InvivoGen provides an untagged and non-glycosylated recombinant human IL-16 in its mature and intracellular form. Its size, purity, and ability to be detected by a specific anti-IL-16 monoclonal antibody were verified (Figures 1 and 2). Importantly, InvivoGen's recombinant hIL-16 is guaranteed free of bacterial contaminations (e.g. lipoproteins and endotoxins) which could interfere with IL-16 specific signaling in TLR2- and TLR4-expressing target cells (Figure 3).

## Detection by SDS-PAGE and WES

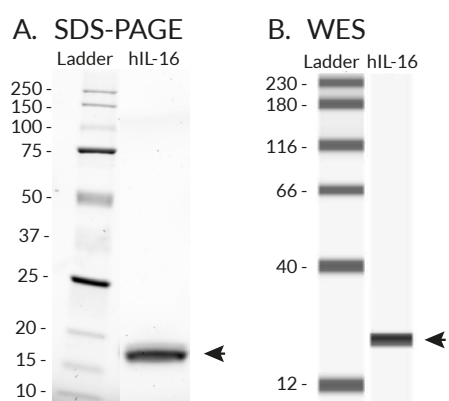


Figure 1: SDS PAGE and WES analysis of the recombinant human (h)IL-16 protein. (A) 0.5  $\mu$ g of hIL-16 was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. A band was detected at ~17 kDa. (B) 4  $\mu$ g of hIL-16 was analyzed by WES using InvivoGen's anti-hIL-16 hIgG1 antibody, followed by an HRP-conjugated anti-human secondary antibody. A band was detected at ~23 kDa.

## Detection by ELISA

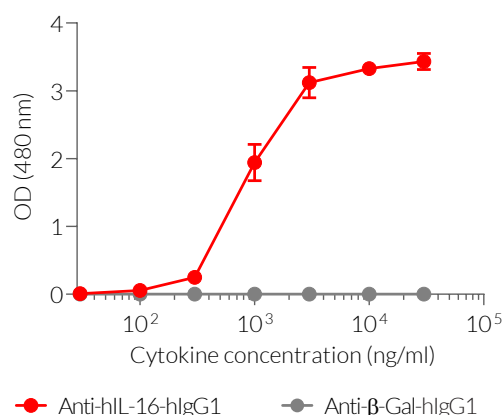


Figure 2: Detection of the recombinant human (h)IL-16 protein by ELISA. Coated ELISA plates with increasing concentrations of hIL-16 were incubated with 2  $\mu$ g/ml of Anti-hIL-16-hIgG1 mAb (red curve) or Anti- $\beta$ -Gal-hIgG1 control antibody (grey curve) for 2h, followed by an HRP-conjugated anti-human secondary antibody (1/1000 dilution). Data are shown as optical density (OD) at 480 nm (mean + SEM).

## Absence of bacterial contamination in recombinant hIL-16

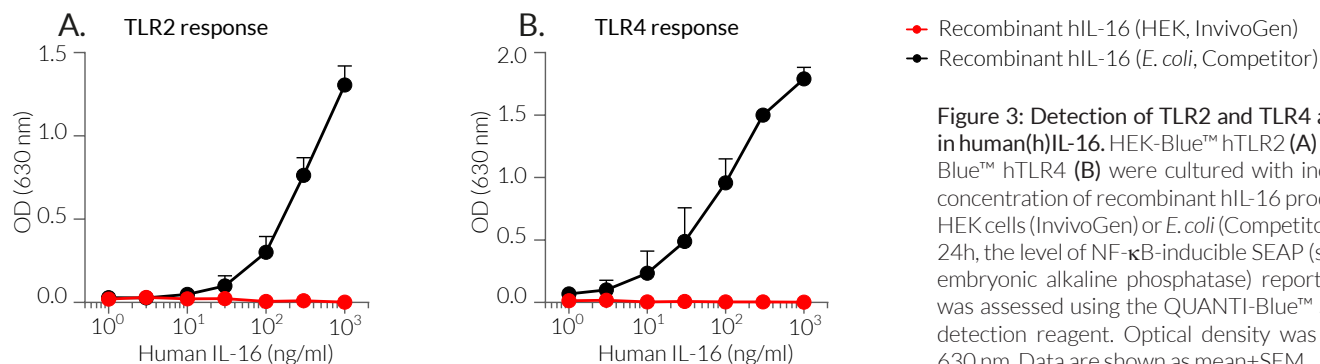


Figure 3: Detection of TLR2 and TLR4 agonists in human(h)IL-16. HEK-Blue™ hTLR2 (A) or HEK-Blue™ hTLR4 (B) were cultured with increasing concentration of recombinant hIL-16 produced in HEK cells (InvivoGen) or *E. coli* (Competitor). After 24h, the level of NF- $\kappa$ B-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene was assessed using the QUANTI-Blue™ Solution detection reagent. Optical density was read at 630 nm. Data are shown as 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)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)