

Validation data for Anti-HER2-hlgG1

<https://www.invivogen.com/anti-her2-0>

For research use only

Version 23J09-NJ

Anti-HER2-hlgG1 has been specifically developed to generate antibody-drug conjugates (ADCs). The antibody specificity for HER2 has been validated using cell surface staining and analysis by flow cytometry (Figure 1). The biological activity of ADCs combining Anti-HER2-hlgG1 and STG-982, a STING agonist, or TL7-887, a TLR7 agonist, has been validated using cellular assays. In a co-culture of HER2⁺ tumor cells (SK-BR-3) and human peripheral blood monocytes (PBMCs), Anti-HER2/STG-982 and Anti-HER2/TL7-887 induce a significantly higher production of CXCL10 and IL-6, respectively, than unconjugated agonists or negative control ADCs (Figures 2A and 3A). Of note, in absence of tumor cells, PBMCs remain either unresponsive to all stimulations (Figure 2B), or respond to higher doses of Anti-HER2/TL7-887 (Figure 3B). This latter observation could be explained by cellular uptake through endocytosis after binding to Fc-receptor and/or to HER2 which is expressed at low levels by some PBMC subpopulations¹.

1. You F. et al., 2008. Low-level expression of HER2 and CK19 in normal peripheral blood mononuclear cells: relevance for detection of circulating tumor cells. *J Hematol Oncol.* 1(2): DOI: 10.1186/1756-8722-1-2.

Validation of Anti-HER2-hlgG1 by flow cytometry

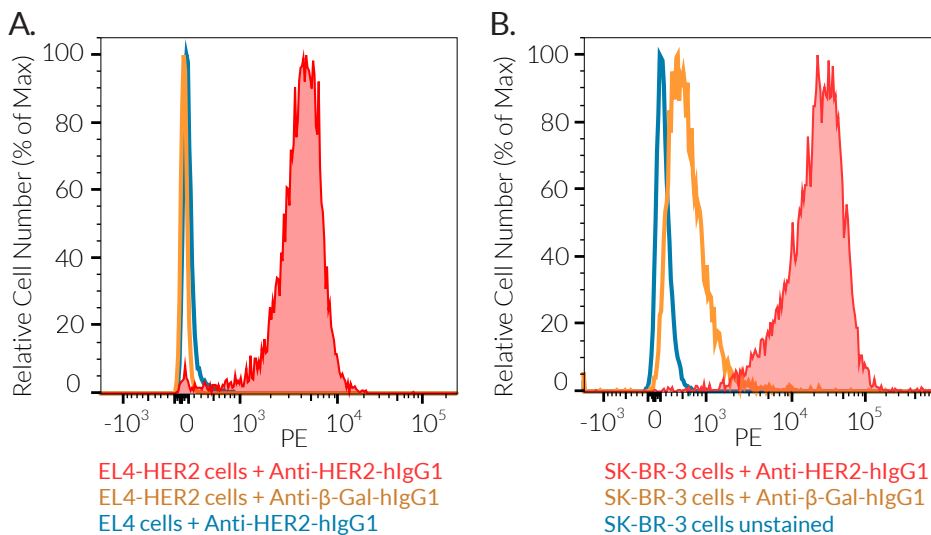


Figure 1: Cell surface staining of HER2 using Anti-HER2-hlgG1 mAb.

~5 × 10⁵ EL4-HER2 (A) or ~3 × 10⁵ SK-BR-3 (B) cells were incubated with 500 ng (A) or 1 μg (B) of Anti-HER2-hlgG1 mAb or an isotype control for 30 min to 1h at 4°C. Cells were then washed and incubated with 250 ng of goat anti-human κ light chain antibody coupled to PE for 45 min to 1h at 4°C. Cell surface staining was analyzed by flow cytometry.

Biological activity of Anti-HER2/STG-982 in co-cultures

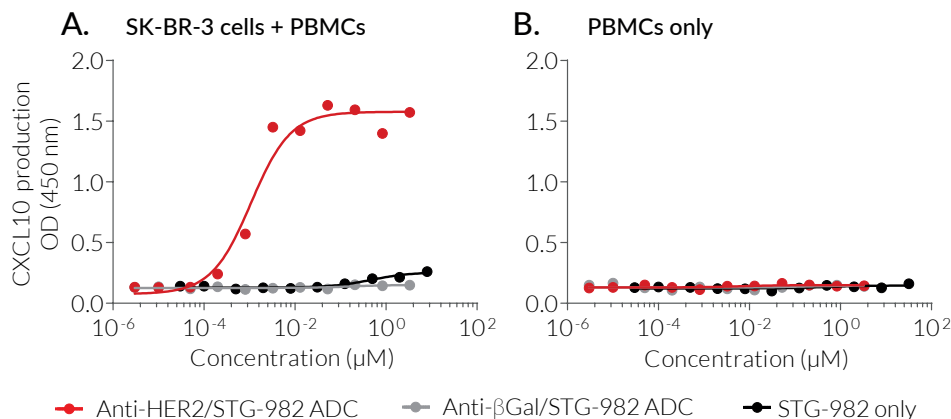


Figure 2: Dose-response of human PBMCs co-cultured with SK-BR-3 tumor cells and Anti-HER2/STG-982 ADC.

1.5 × 10⁵ human PBMCs and 5 × 10⁴ SK-BR-3 tumor cells (A) or 1.5 × 10⁵ human PBMCs only (B) were incubated with increasing concentrations of Anti-HER2/STG-982 ADC (DAR ~5), Anti-β-Gal/STG-982 ADC (DAR ~5), or STG-982 only. After overnight incubation, the STING-mediated response was assessed by measuring the production of CXCL10 in PBMC and SK-BR-3 co-culture supernatants, using an ELISA. The optical density (OD) at 450 nm is shown.

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

Biological activity of Anti-HER2/TL7-887 in co-cultures

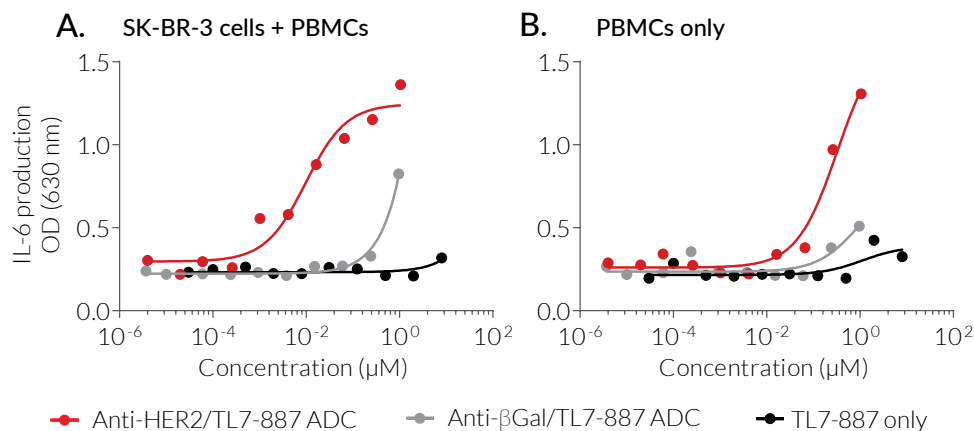


Figure 3: Dose-response of human PBMCs co-cultured with SK-BR-3 tumor cells and Anti-HER2/TL7-887 ADC.

1.5×10^5 human PBMCs and 5×10^4 SK-BR-3 tumor cells (A) or 1.5×10^5 human PBMCs only (B) were incubated with increasing concentrations of Anti-HER2/TL7-887 ADC (DAR ~6), Anti- β -Gal/TL7-887 ADC (DAR ~6), or TL7-887 only. After overnight incubation, the TLR7-mediated response was determined using HEK-Blue™ IL-6 reporter cells. Briefly, the levels of IL-6 production in PBMC and SK-BR-3 co-culture supernatants were assessed by measuring the SEAP activity of HEK-Blue™ IL-6 reporter cells, using QUANTI-Blue™, a SEAP detection reagent. The optical density (OD) at 630 nm is shown.

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