

Validation data for Anti-TROP2-hlgG1

<https://www.invivogen.com/anti-trop2>

For research use only

Version 23J09-NJ

Anti-TROP2-hlgG1 has been specifically developed to generate antibody-drug conjugates (ADCs). The antibody specificity for TROP2 has been validated using flow cytometry and ELISA (Figure 1). The biological activity of ADCs combining Anti-TROP2-hlgG1 and STG-982, a STING agonist, or TL7-887, a TLR7 agonist, has been validated using cellular assays. In a co-culture of TROP2⁺ tumor cells (BxPC-3) and human peripheral blood monocytes (PBMCs), Anti-TROP2/STG-982 and Anti-TROP2/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- β -gal/TL7-887 control ADC (Figure 3B), which could be explained by cellular uptake through endocytosis/pinocytosis.

Validation of Anti-TROP2-hlgG1 by flow cytometry and ELISA

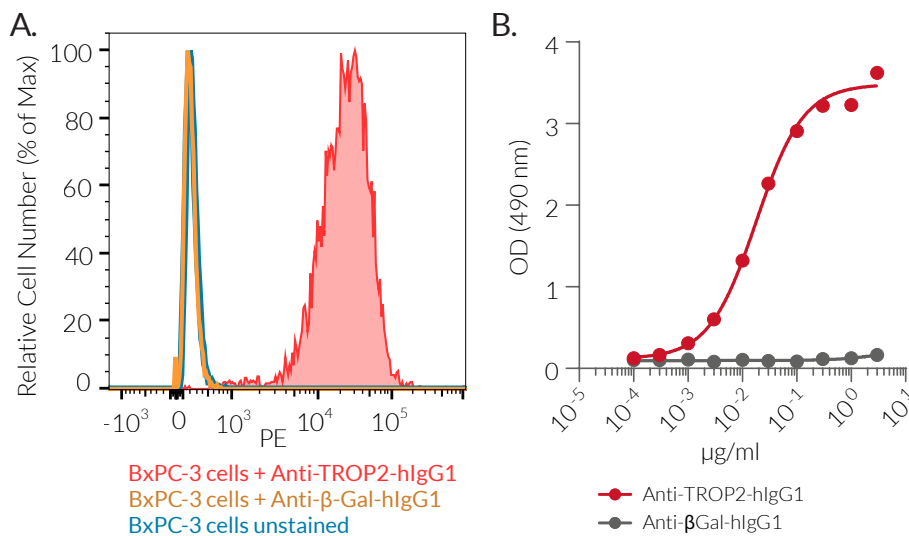


Figure 1: Anti-TROP2-hlgG1 mAb binding to cell-surface TROP2 or plate-bound TROP2.

(A) $\sim 5 \times 10^5$ BxPC-3 tumor cells were incubated with 1 μ g of Anti-TROP2-hlgG1 mAb or an isotype control for 45 min at 4°C. Cells were then washed and incubated with 250 ng of goat anti-human κ light chain antibody coupled to PE for 1h at 4°C. Cell surface staining was analyzed by flow cytometry.

(B) TROP2-Fc fusion protein (1 μ g/ml) was coated on ELISA plates overnight. A 3-fold serial dilution of Anti-TROP2-hlgG1 (red curve) or of Anti- β -Gal-hlgG1 control mAb (grey curve) was realized for the capture step. An HRP-labelled anti-human κ light chain antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 490 nm.

Biological activity of Anti-TROP2/STG-982 in co-cultures

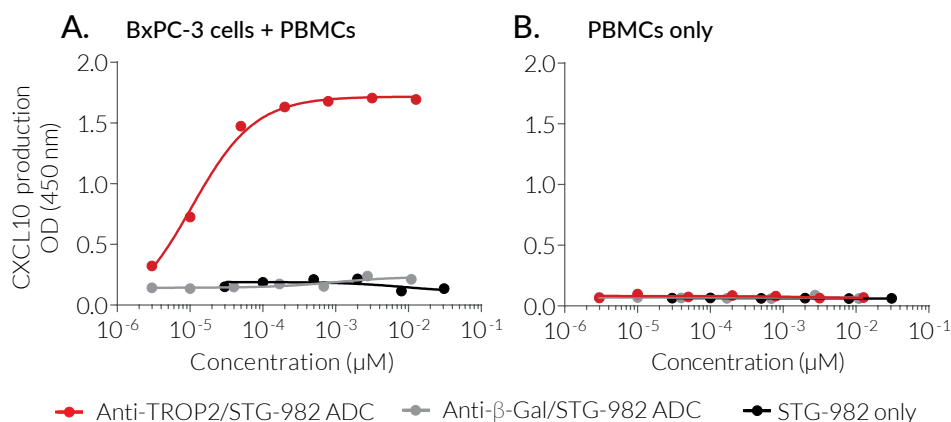


Figure 2: Dose-response of human PBMCs co-cultured with BxPC-3 tumor cells and Anti-TROP2/STG-982 ADC.

1.5×10^5 human PBMCs and 5×10^4 BxPC-3 tumor cells (A) or 1.5×10^5 human PBMCs only (B) were incubated with increasing concentrations of Anti-TROP2/STG-982 ADC (DAR ~ 4), Anti- β -Gal/STG-982 ADC (DAR ~ 4), or STG-982 only. After overnight incubation, the STING-mediated response was assessed by measuring the production of CXCL10 in PBMC and BxPC-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-TROP2/TL7-887 in co-cultures

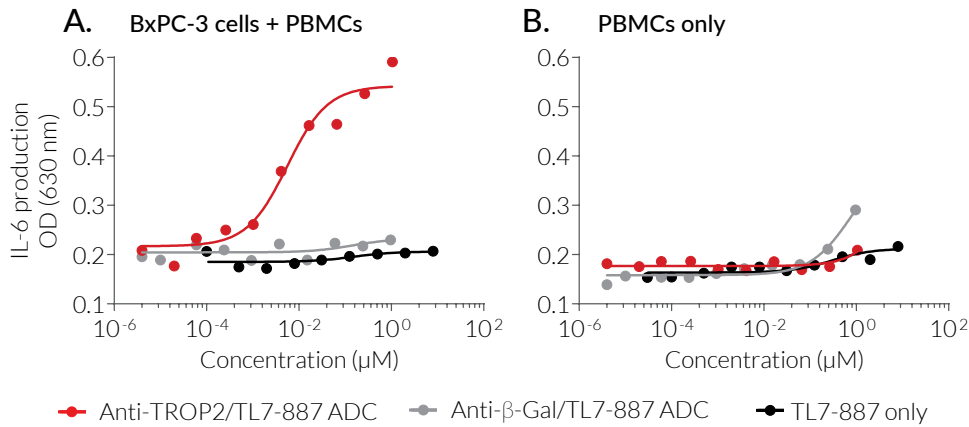


Figure 3: Dose-response of human PBMCs co-cultured with BxPC-3 tumor cells and Anti-TROP2/TL7-887 ADC.

1.5 × 10⁵ human PBMCs and 5 × 10⁴ BxPC-3 tumor cells **(A)** or 1.5 × 10⁵ human PBMCs only **(B)** were incubated with increasing concentrations of Anti-TROP2/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 BxPC-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