

Validation data for Anti-β-Gal-hlgG1*

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

For research use only

Version 23J12-NJ

Anti-β-Gal-hlgG1* has been specifically developed for the generation of antibody-drug conjugates (ADCs). This antibody can be used to generate control ADCs which do not recognize any antigen other than β-galactosidase. The antibody specificity for β-galactosidase has been validated using ELISA (Figure 1). The absence of non-specific binding and biological activity of Anti-β-Gal-hlgG1* has been validated using flow cytometry (Figure 2) and antibody-drug conjugate (ADC)-based cellular assays (Figure 3), respectively.

Validation of Anti-β-Gal-hlgG1* by ELISA

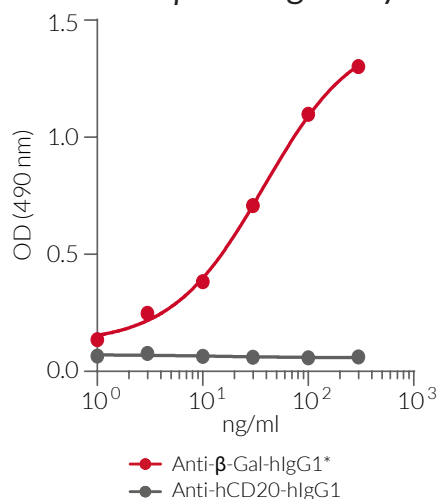


Figure 1: Anti-β-hlgG1* mAb binding to plate-bound β-galactosidase. β-galactosidase protein (2 μg/ml) was coated on ELISA plates overnight. A 3-fold serial dilution of Anti-β-Gal-hlgG1* (red curve) or of Anti-hCD20-hlgG1 control mAb (grey curve) was realized for the capture step. An HRP-labelled anti-human IgG1 (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 490 nm.

Validation of Anti-β-Gal-hlgG1* by flow cytometry

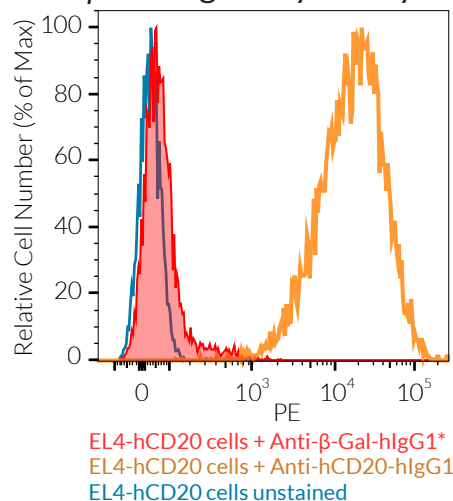


Figure 2: Absence of Anti-β-hlgG1* mAb binding to EL4-hCD20 cells. ~5 × 10⁵ EL4 cells stably expressing cell surface human CD20 antigen were incubated with 2 μg of Anti-β-Gal-hlgG1* mAb or Anti-hCD20-hlgG1 control mAb for 45 min at 4°C. Cells were then washed and incubated with 250 ng of goat anti-human κ light chain antibody coupled to PE for 1h at 4°C. Cell surface staining was analyzed by flow cytometry.

Absence of biological activity of Anti-β-Gal/STG-982 in co-cultures

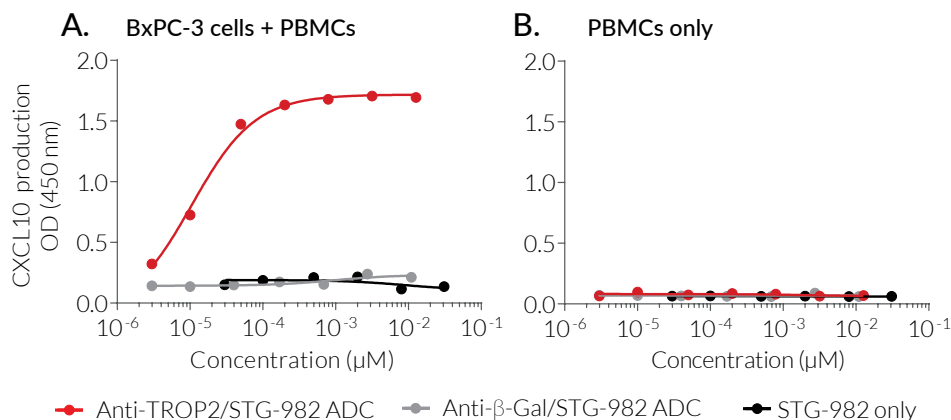


Figure 2: Dose-response of human PBMCs co-cultured with BxPC-3 tumor cells and Anti-TROP2/STG-982 or Anti-β-Gal/STG982 ADC.

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/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