

Validation data for Raji-HER2 Cells

<https://www.invivogen.com/raji-her2>

For research use only

Version 21B17-ED

Raji-HER2 cells were developed from the Raji cell line to overexpress the human HER2 gene. Raji-HER2 cells were designed as target cells in InvivoGen's antibody-dependent cellular cytotoxicity (ADCC) assay using clinically-relevant anti-HER2 monoclonal antibodies (mAbs). HER2 expression by Raji-HER2 cells has been verified by flow-cytometry (Figure 1), and induction of ADCC has been validated using InvivoGen's collection of anti-HER2 antibody isotypes and Jurkat-Lucia™ NFAT-CD16 reporter cells (Figure 2). The level of ADCC induction is measured by an NFAT-dependent Lucia luciferase reporter protein.

Validation of HER2 expression by flow cytometry

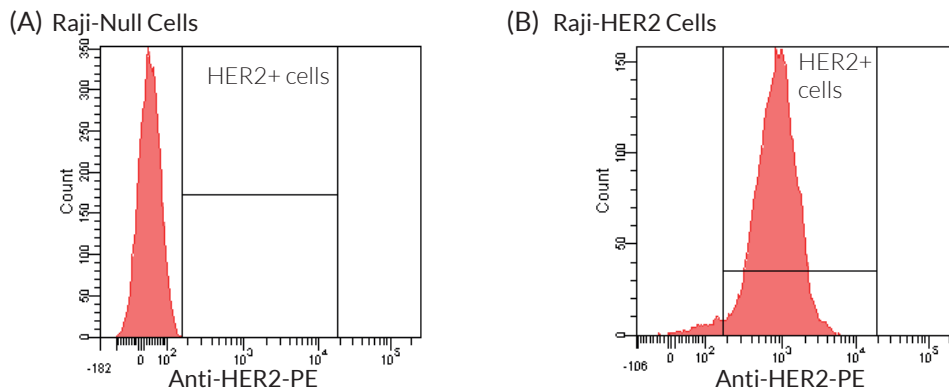


Figure 1: Validation of the expression of HER2 by Raji-HER2 cells. Raji-Null (A) and Raji-HER2 (B) cells were incubated with a PE-conjugated Anti-HER2 mAb for 30 minutes. The binding affinity was then measured using flow cytometry.

ADCC assay using various anti-HER2 antibody isotypes and Raji-HER2 target cells

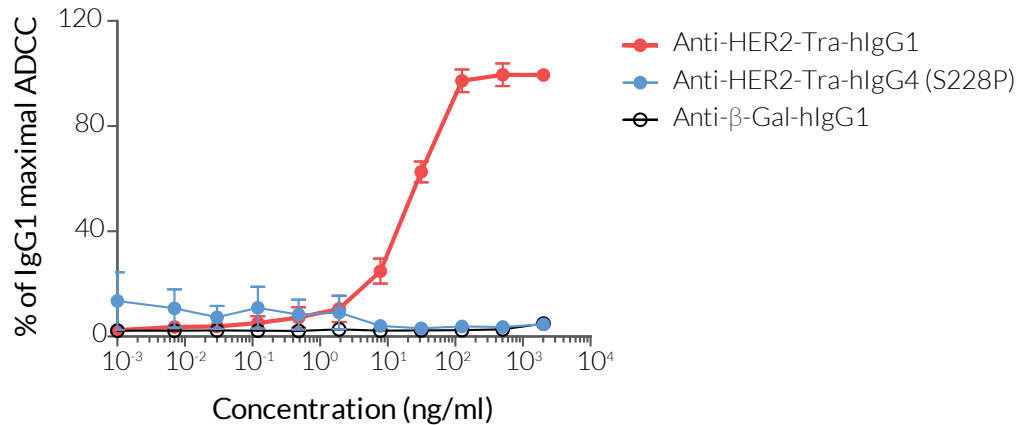


Figure 2: Comparison of ADCC potency for native and engineered anti-HER2 antibody isotypes. Raji-HER2 cells were incubated with gradient concentrations of Anti-HER2-Tra mAbs, which feature the Trastuzumab variable region, or an Anti-β-galactosidase (β-Gal) mAb for 1 hour. Jurkat-Lucia™ NFAT-CD16 effector cells were then co-incubated with target cells for 6 hours. NFAT activation, reflecting the induced ADCC response, was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Percentages of the maximal response normalized to the IgG1 isotype are 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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com