

# Validation data for CTLA4/CD80 Bio-IC™

[invivogen.com/immune-checkpoint-hctla4-hcd80-bioassay](http://invivogen.com/immune-checkpoint-hctla4-hcd80-bioassay)

For research use only

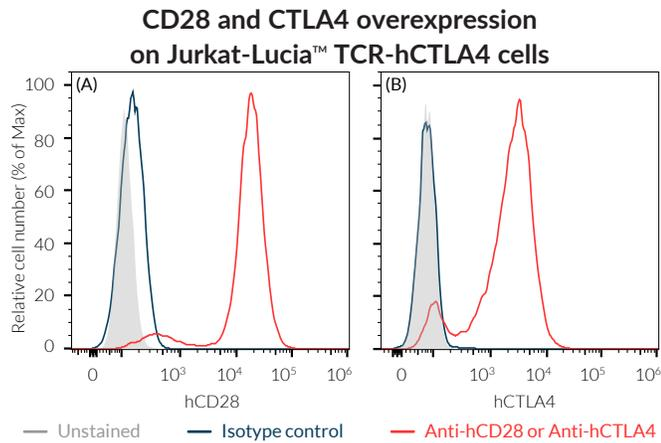
Version 25F12-NJ

CTLA4/CD80 Bio-IC™ is a bioluminescent cell-based assay designed for the screening of novel inhibitors of the CTLA4/CD80 immune checkpoint (IC) axis, such as monoclonal antibodies (mAbs), Fc-fusion proteins, or small molecule-based inhibitors. The assay relies on the co-culture of two cell lines:

**Jurkat-Lucia™ TCR-hCTLA4 effector cells** stably express a specific [HLA::peptide]-restricted TCR and an NFAT-inducible Lucia luciferase reporter gene. In addition, these cells overexpress the human (h)CD28 (cluster of differentiation 28) stimulatory receptor and the hCTLA4 (cytotoxic T lymphocyte-associated protein 4) inhibitory receptor (**Figure 1**).

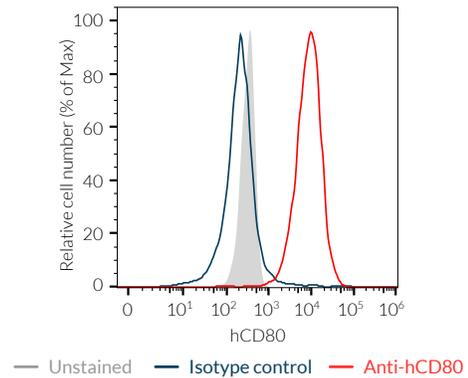
**Raji-APC-Null cells** stably express the specific [HLA::peptide]. They express endogenous hCD80, the inhibitory ligand for CTLA4 (**Figure 2**).

This cellular assay has been functionally tested using the biosimilar Anti-hCTLA4 Ipilimumab. Upon addition of the blocking mAb, the CTLA4/CD80 inhibitory interaction is disrupted and Jurkat-Lucia™ TCR-hCTLA4 effector cells express Lucia luciferase in a dose-dependent manner (**Figure 3**).



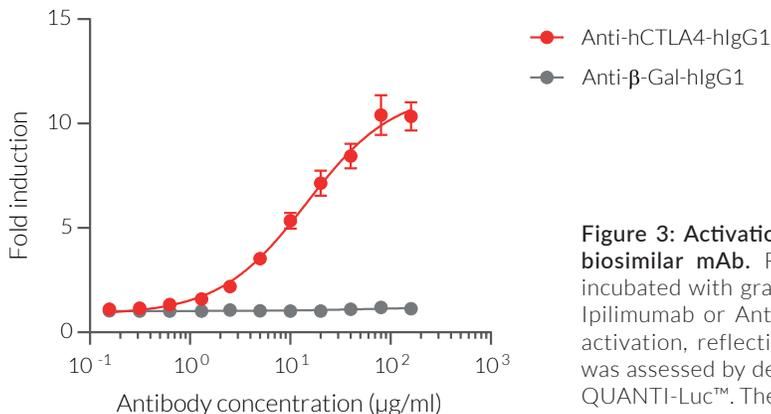
**Figure 1: Validation of human CD28 and CTLA4 overexpression by Jurkat-Lucia™ TCR-hCTLA4 cells.** Jurkat-Lucia™ TCR-hCTLA4 cells were incubated with a PE-conjugated Anti-hCD28 (A) or APC-conjugated Anti-hCTLA4 (B) mAb for 30 minutes. The binding affinity was then measured using flow cytometry.

## Endogenous CD80 expression on Raji-APC-Null cells



**Figure 2: Validation of endogenous human CD80 expression by Raji-APC-Null cells.** Raji-APC-Null cells were incubated with a PE-conjugated Anti-hCD80 mAb for 30 minutes. The binding affinity was then measured using flow cytometry.

## Disruption of CTLA4/CD80 inhibitory interaction using Anti-hCTLA4 biosimilar antibody



**Figure 3: Activation of Jurkat-Lucia™ TCR-hCTLA4 cells using Anti-hCTLA4 biosimilar mAb.** Raji-APC-Null and Jurkat-Lucia™ TCR-hCTLA4 cells were incubated with gradient concentrations of the biosimilar Anti-hCTLA4-hIgG1 Ipilimumab or Anti-β-Gal-hIgG1, as a negative control, for 24 hours. NFAT activation, reflecting the disruption of CTLA4/CD80 inhibitory interaction, was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. The fold increase over non induced cells (no mAbs) 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](mailto:info@invivogen.com)

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