

Validation data for Jurkat-Lucia™ NFAT-CD32 Cells

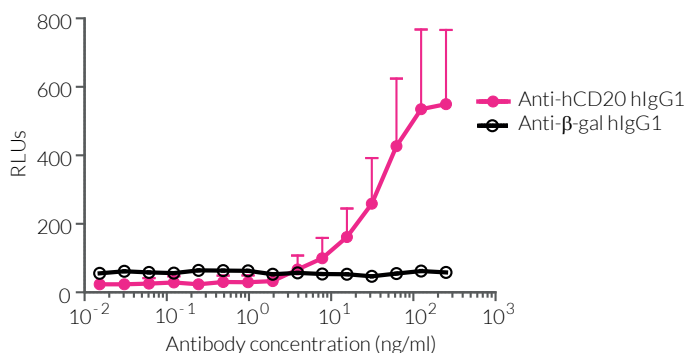
<https://www.invivogen.com/jurkat-lucia-nfat-adcc-adcp-cells>

For research use only

Version 21C16-NJ

Jurkat-Lucia™ NFAT-CD32 cells were engineered from the human T-lymphocyte Jurkat cell line and designed as reporter cells for antibody-dependent cell-mediated phagocytosis (ADCP). Nuclear translocation of NFAT (nuclear factor of activated T-cells), a transcription factor naturally expressed by Jurkat cells, is an early signaling event in ADCP induction. Jurkat-Lucia™ NFAT-CD32 cells stably express the cell surface Fc receptor CD32A (FcγRIIA; H1611 allotype) and the Lucia luciferase reporter gene under the control of an ISG54 minimal promoter fused to six NFAT response elements. Human CD32A expression by Jurkat-Lucia™ NFAT-CD32 cells has been verified by flow-cytometry. These cells have been functionally tested with various target cells and specific monoclonal antibody (mAb) isotype combinations.

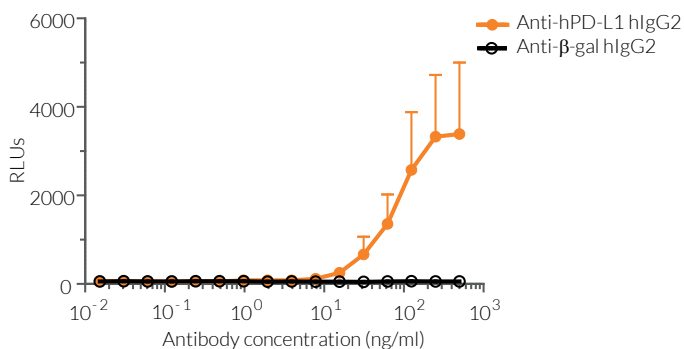
Jurkat-Lucia™ NFAT-CD32 cell responses to ADCP induction with Anti-hCD20 hIgG1 and Raji-Null cells



ADCP potency of Anti-hCD20 hIgG1:

Raji-Null cells (expressing hCD20) were incubated with gradient concentrations of Anti-hCD20 hIgG1 (featuring the variable region of Rituximab) or Anti-β-galactosidase (β-gal) hIgG1 for 1 hour. Jurkat-Lucia™ NFAT-CD32 effector cells were then co-incubated with target cells for 6 hours. NFAT activation, reflecting the induced ADCP response, was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Relative light units (RLUs) are shown.

Jurkat-Lucia™ NFAT-CD32 cell responses to ADCP induction with Anti-hPD-L1 hIgG2 and Raji-hPD-L1 cells



ADCP potency of Anti-hPD-L1 hIgG2:

Raji-hPD-L1 cells were incubated with gradient concentrations of Anti-hPD-L1 hIgG2 or Anti-β-galactosidase (β-gal) hIgG2 for 1 hour. Jurkat-Lucia™ NFAT-CD32 effector cells were then co-incubated with target cells for 6 hours. NFAT activation, reflecting the induced ADCP response, was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Relative light units (RLUs) 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



Any questions about our cell lines?
Visit our FAQ page.

 **InvivoGen**
www.invivogen.com