

Validation data for Jurkat-Lucia™ NFAT Cells

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

For research use only

Version 22C03-NJ

Jurkat-Lucia™ NFAT cells were designed from a clone of the human T lymphocyte Jurkat cell line deficient for CD28 expression (Figure 1). These cells stably express the Lucia luciferase reporter gene under the control of an ISG54 minimal promoter fused to six NFAT response elements. The activation of NFAT has been confirmed following the induction of Jurkat-Lucia™ NFAT cells with concanavalin A, phytohaemagglutinin P, or PMA/ionomycin (Figure 2).

CD3 and CD28 expression on Jurkat-Lucia™ NFAT cells

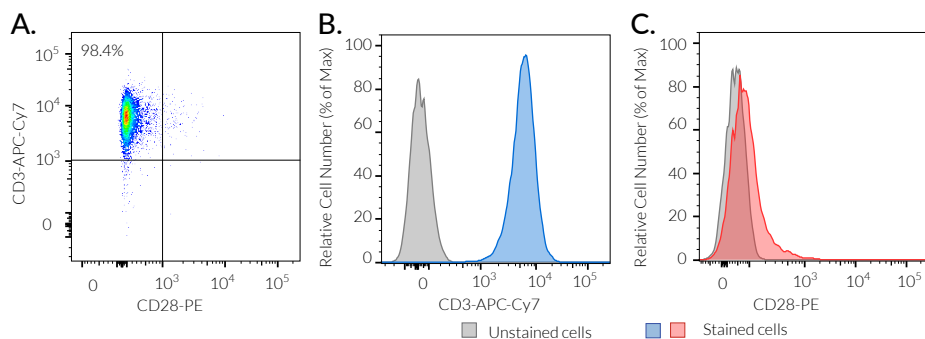


Figure 1: Validation of human CD3 expression and absence of CD28 expression by Jurkat-Lucia™ NFAT cells. Jurkat-Lucia™ NFAT cells were incubated with a cocktail of APC-Cy7-conjugated anti-hCD3 and PE-conjugated anti-hCD28 antibodies for 30 minutes. The binding affinity was then measured using flow-cytometry. (A) CD3 and CD28 co-staining. (B) CD3 and (C) CD28 expression compared to unstained cells.

Jurkat-Lucia™ NFAT cell responses

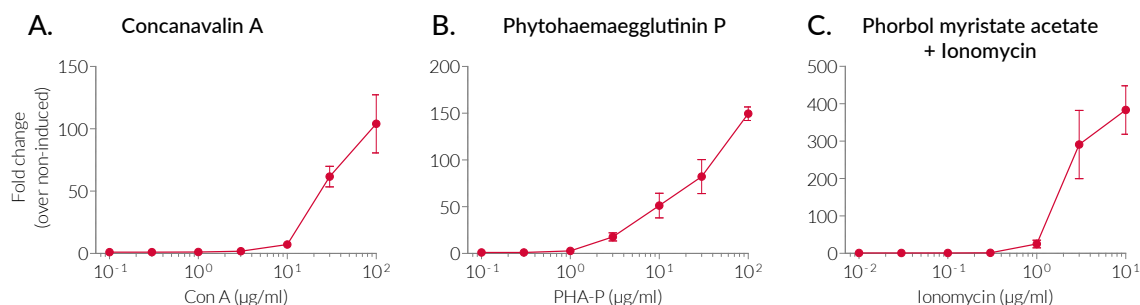


Figure 2: Validation of NFAT activation in Jurkat-Lucia™ NFAT cells.

Jurkat-Lucia™ NFAT cells were incubated with increasing concentrations of concanavalin A (Con A) (A), phytohaemagglutinin P (PHA-P) (B), or Ionomycin with 50 ng/ml phorbol myristate acetate (PMA) (C) for 24 hours. NFAT activation was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Fold change over non-induced cells 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