

Validation data for Jurkat-Dual™ Cells

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

For research use only

Version 21L22-MM

Jurkat-Dual™ cells were derived from the human T lymphocyte-based Jurkat cell line by the stable integration of two inducible reporter constructs: Lucia luciferase and SEAP (secreted embryonic alkaline phosphatase). As a result, they allow the simultaneous study of the NF- κ B pathway, by assessing the activity of Lucia luciferase, and the IRF (interferon regulatory factor) pathway, by monitoring the activity of SEAP.

Jurkat-Dual™ cells induce the activation of NF- κ B and produce Lucia luciferase in response to TNF- α and T-lymphocyte mitogens, such as phytohemagglutinin (PHA) and concanavalin A (ConA). Additionally, they trigger the IRF pathway and produce SEAP upon stimulation with type I interferons (IFNs) or poly(I:C) (Figure 1).

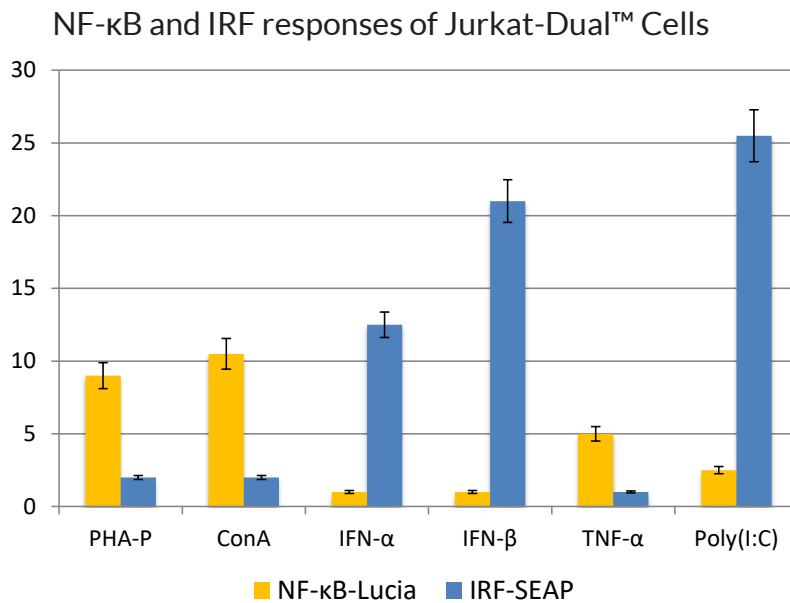


Figure 1. NF- κ B/IRF dual response of Jurkat-Dual™ cells to various stimuli. Cells were incubated with 20 μ g/ml of the mitogens phytohemagglutinin-P (PHA-P, the protein form of PHA) or concanavalin A (ConA), 300 IU/ml IFN- α , 300 IU/ml IFN- β , 100 ng/ml TNF- α or 10 μ g/ml poly(I:C). After 24h incubation, NF- κ B and IRF activation was assessed by measuring the levels of SEAP and Lucia luciferase in the supernatant using QUANTI-Luc™ or QUANTI-Blue™ Solution, respectively.

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