Validation data for THP1-Dual[™] KO-cGAS cells

https://www.invivogen.com/thp1-dual-kocgas

For research use only

Version 19K16-MM

THP1-DualTM KO-cGAS cells were generated from THP1-DualTM cells by stable knockout of the cGAS gene which has been confirmed by Western blot (figure 1). These cells derive from the human monocytic cell line THP-1, which is often used for the study of DNA-sensing pathways. THP1-DualTM and THP1-DualTM KO-cGAS cells can be used to study cGAS signaling. Both cell lines express two inducible reporter constructs that enable the simultaneous study of the NF- κ B pathway, by monitoring the activity of SEAP (secreted embryonic alkaline phosphatase), and the IRF (interferon regulatory factor) pathway, by assessing the activity of the secreted Lucia luciferase. The IRF induction response of these cells to different ligands has been assessed (figure 2). RAW-LuciaTM ISG-KO-cGAS cells respond to interferons (e.g. IFN- α and IFN- β) and cyclic dinucelotides (e.g. 2'3'-cGAMP). However, as expected, they respond very poorly to transfected DNA, such as VACV70/LyoVecTM and poly(dA:dT)/LyoVecTM. The NF- κ B response of THP1-DualTM KO-cGAS cells is unaffected by the knockout of the cGAS gene (figure 3).





NF-κB Induction (SEAP reporter)



Figure 2: Stimulation of THP1-DualTM (parental cell line) and THP1-DualTM KO-cGAS cells with VACV70/LyoVecTM (1 µg/ml), poly(dA:dT)/LyoVecTM (100 ng/ml), and 2'3'-cGAMP (1 µg/ml). Human IFN- α (1 x 10⁴ U/ml) and IFN- β (1 x 10⁴ U/ml) serve as positive controls. Human TNF- α (10 ng/ml) has been included as a negative control. After a 24h incubation, IRF activation was determined by measuring the relative light units (RLUs) in a luminometer using QUANTI-LucTM, a Lucia luciferase detection reagent. The IRF induction of each ligand is expressed relative to that of hIFN- β at 1 x 10⁴ U/ml (taken as 100%). Figure 3: THP1-DualTM and THP1-DuaTM KO-cGAS cells were incubated with TNF- α (0.1 ng/ml), Pam3CSK4 (0.1 ng/ml; TLR1/2 ligand), LPS-EB Ultrapure (100 ng/ml; TLR4 ligand), VACV70/LyoVecTM (1 µg/ml), poly(dA:dT)/LyoVecTM (100 ng/ml), 2'3'-cGAMP (3 µg/ml) and hIFN- α (1 x 10⁴ U/ml). After a 24h incubation, NF- κ B activation was determined using QUANTI-BlueTM, a SEAP detection reagent, and by reading the optical density (OD) at 655 nm.

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

